

---

## PMA & PMAxx™ Selected References

[Click on the blue link to see journal article](#)

### Contents:

#### Viability PCR Reviews

p. 1

#### Viability PCR Dyes

- PMAxx™
- PMA

p. 2

There are hundreds of journal articles using PMA in viability PCR. Publications using PMA in select applications and cell types have been selected below.

#### Related Products

- PMA-Lite™
- PMA Enhancer

p. 2

p. 3

#### Sample Types

- Bacteria

The use of PMA for viability PCR in bacteria has been published in hundreds of journal articles. See select articles below for certain applications or techniques.

- Yeast and Fungi
- Biofilms
- Eukaryotes
- Viruses
- Archaea

p. 3

p. 4

p. 6

p. 6

p. 9

#### Applications

- Food Science
- Probiotics
- Environmental Testing
- Filtered Samples

p. 9

p. 13

p. 14

p. 17

#### Techniques

- Sequencing
- Loop-mediated Isothermal Amplification (LAMP)

p. 17

p. 20

### Viability PCR Reviews

Elizaquivel, P., Aznar, R., and Sanchez, G. (2013). [Recent developments in the use of viability dyes and quantitative PCR in the food microbiology field](#). J Appl Microbiol 116, 1-13. 10.1111/jam.12365

Emerson, J. B., Adams, R. I., Roman, C. M. B., Brooks, B., Coil, D. A., Dahlhausen, K., Ganz, H. H., Hartmann, E. M., Hsu, T., Justice, N. B., et al. (2017). [Schrodinger's microbes: Tools for distinguishing the living from the dead in microbial ecosystems](#). Microbiome 5, 86. 10.1186/s40168-017-0285-3

Fittipaldi, M., Nocker, A., and Codony, F. (2012). [Progress in understanding preferential detection of live cells using viability dyes in combination with DNA amplification](#). J Microbiol Methods 91, 276-289. 10.1016/j.mimet.2012.08.007

Greening, G. (2013). [Foodborne viruses: a focus on challenges associated with detection methods](#). Microbiology Australia DOI: 10.1071/MA13022. 10.1071/MA13022

Li, Y., Yang, L., Fu, J., Yan, M., Chen, D., and Zhang, L. (2017). [The novel loop-mediated isothermal amplification based confirmation methodology on the bacteria in Viable but Non-Culturable \(VBNC\) state](#). *Microb Pathog* 111, 280-284. S0882-4010(17)30999-3

van Frankenhuyzen, J. K., Trevors, J. T., Lee, H., Flemming, C. A., and Habash, M. B. (2011). [Molecular pathogen detection in biosolids with a focus on quantitative PCR using propidium monoazide for viable cell enumeration](#). *J Microbiol Methods* 87, 263-272. 10.1016/j.mimet.2011.09.007

#### **PMAxx™**

Emerson, J. B., Adams, R. I., Roman, C. M. B., Brooks, B., Coil, D. A., Dahlhausen, K., Ganz, H. H., Hartmann, E. M., Hsu, T., Justice, N. B., et al. (2017). [Schrodinger's microbes: Tools for distinguishing the living from the dead in microbial ecosystems](#). *Microbiome* 5, 86. 10.1186/s40168-017-0285-3

Fraisse, A., Niveau, F., Hennechart-Collette, C., Coudray-Meunier, C., Martin-Latil, S., and Perelle, S. (2018). [Discrimination of infectious and heat-treated norovirus by combining platinum compounds and real-time RT-PCR](#). *Int J Food Microbiol* 269, 64-74. S0168-1605(18)30023-0

Garcia-Fontana, C., Narvaez-Reinaldo, J. J., Castillo, F., Gonzalez-Lopez, J., Luque, I., and Manzanera, M. (2016). [A New Physiological Role for the DNA Molecule as a Protector against Drying Stress in Desiccation-Tolerant Microorganisms](#). *Front Microbiol* 7, 2066. 10.3389/fmicb.2016.02066

Randazzo, W., Lopez-Galvez, F., Allende, A., Aznar, R., and Sanchez, G. (2016). [Evaluation of viability PCR performance for assessing norovirus infectivity in fresh-cut vegetables and irrigation water](#). *Int J Food Microbiol* 229, 1-6. S0168-1605(16)30169-6

Randazzo, W., Piqueras, J., Rodriguez-Diaz, J., Aznar, R., and Sanchez, G. (2018). [Improving efficiency of viability-qPCR for selective detection of infectious HAV in food and water samples](#). *J Appl Microbiol* 124, 958-964. 10.1111/jam.13519

#### **PMA-Lite™**

Burnet, J. B., Faraj, T., Cauchie, H. M., Joaquim-Justo, C., Servais, P., Prevost, M., and Dorner, S. M. (2017). [How does the cladoceran \*Daphnia pulex\* affect the fate of \*Escherichia coli\* in water?](#) *PLoS One* 12, e0171705. 10.1371/journal.pone.0171705

Eramo, A., Medina, W. M., and Fahrenfeld, N. L. (2017). [Peracetic acid disinfection kinetics for combined sewer overflows: indicator organisms, antibiotic resistance genes, and microbial community](#). *Environ Sci (Camb)* 3, 1061-1072. 10.1039/C7EW00184C

Oh, E., McMullen, L., and Jeon, B. (2015). [Impact of oxidative stress defense on bacterial survival and morphological change in \*Campylobacter jejuni\* under aerobic conditions](#). *Front Microbiol* 6, 295. 10.3389/fmicb.2015.00295

Overney, A., Jacques-Andre-Coquin, J., Ng, P., Carpentier, B., Guillier, L., and Firmesse, O. (2017). [Impact of environmental factors on the culturability and viability of \*Listeria monocytogenes\* under conditions encountered in food processing plants](#). *Int J Food Microbiol* 244, 74-81. 10.1016/j.ijfoodmicro.2016.12.012

Yang, Y., Cheng, D., Li, Y., Yu, L., Gin, K. Y., Chen, J. P., and Reinhard, M. (2017). [Effects of monochloramine and hydrogen peroxide on the bacterial community shifts in biologically treated wastewater](#). *Chemosphere* 189, 399-406. 10.1016/j.chemosphere.2017.09.087

Xiao, L., Zhang, Z., Sun, X., Pan, Y., and Zhao, Y. (2015). [Development of a quantitative real-time PCR assay for viable \*Salmonella\* spp. without enrichment](#). *Food Control*. 10.1016/j.foodcont.2015.03.050

Zhang, Z., Liu, H., Lou, Y., Xiao, L., Liao, C., Malakar, P. K., Pan, Y., and Zhao, Y. (2015). [Quantifying viable \*Vibrio parahaemolyticus\* and \*Listeria monocytogenes\* simultaneously in raw shrimp](#). *Appl Microbiol Biotechnol* 99, 6451-6462. 10.1007/s00253-015-6715-x

Zheng, Q., Miks-Krajnik, M., D'Souza, C., Yang, Y., Heo, D. J., Kim, S. K., Lee, S. C., and Yuk, H. G. (2015). [Growth of healthy and sanitizer-injured \*Salmonella\* cells on mung bean sprouts in different commercial enrichment broths](#). *Food Microbiol* 52, 159-168. 10.1016/j.fm.2015.07.013

### **PMA Enhancer**

Youn, S. Y., Jeong, O. M., Choi, B. K., Jung, S. C., and Kang, M. S. (2016). [Application of loop-mediated isothermal amplification with propidium monoazide treatment to detect live \*Salmonella\* in chicken carcasses](#). *Poult Sci*. 10.3382/ps/pew341

Duarte-Guevara, P., Duarte-Guevara, C., Ornob, A., and Bashir, R. (2016). [On-chip PMA labeling of foodborne pathogenic bacteria for viable qPCR and qLAMP detection](#). *Microfluid Nanofluid*. 10.1007/s1040

### **Yeast and Fungi**

Agusti, G., Fittipaldi, M., Morato, J., and Codony, F. (2013). [Viable quantitative PCR for assessing the response of \*Candida albicans\* to antifungal treatment](#). *Appl Microbiol Biotechnol* 97, 341-349. 10.1007/s00253-012-4524-z

Andorra, I., Esteve-Zarzoso, B., Guillamon, J. M., and Mas, A. (2010). [Determination of viable wine yeast using DNA binding dyes and quantitative PCR](#). *Int J Food Microbiol* 144, 257-262. 10.1016/j.ijfoodmicro.2010.10.003

Blachowicz, A., Mayer, T., Bashir, M., Pieber, T. R., De Leon, P., and Venkateswaran, K. (2017). [Human presence impacts fungal diversity of inflated lunar/Mars analog habitat](#). *Microbiome* 5, 62. 10.1186/s40168-017-0280-8

Checinska, A., Probst, A. J., Vaishampayan, P., White, J. R., Kumar, D., Stepanov, V. G., Fox, G. E., Nilsson, H. R., Pierson, D. L., Perry, J., et al. (2015). [Microbiomes of the dust particles collected from the International Space Station and Spacecraft Assembly Facilities](#). *Microbiome* 3, 50. 10.1186/s40168-015-0116-3

Crespo-Sempere, A., Estiarte, N., Marin, S., Sanchis, V., and Ramos, A. J. (2013). [Propidium monoazide combined with real-time quantitative PCR to quantify viable \*Alternaria\* spp. contamination in tomato products](#). *Int J Food Microbiol* 165, 214-220. 10.1016/j.ijfoodmicro.2013.05.017

Luo, Y., Bolt, H. L., Eggimann, G. A., McAuley, D. F., McMullan, R., Curran, T., Zhou, M., Jahoda, P. C., Cobb, S. L., and Lundy, F. T. (2016). [Peptoid Efficacy against Polymicrobial Biofilms Determined by Using Propidium Monoazide-Modified Quantitative PCR](#). *Chembiochem*. 10.1002/cbic.201600381

Medina, E., Perez-Diaz, I. M., Breidt, F., Hayes, J., Franco, W., Butz, N., and Azcarate-Peril, M. A. (2016). [Bacterial Ecology of Fermented Cucumber Rising pH Spoilage as Determined by Nonculture-Based Methods](#). *J Food Sci* 81, M121-129. 10.1111/1750-3841.13158

Nguyen, L. D., Deschaght, P., Merlin, S., Loywick, A., Audebert, C., Van Daele, S., Viscogliosi, E., Vaneechoutte, M., and Delhaes, L. (2016). [Effects of Propidium Monoazide \(PMA\) Treatment on Mycobiome and Bacteriome Analysis of Cystic Fibrosis Airways during Exacerbation](#). *PLoS One* 11, e0168860. 10.1371/journal.pone.0168860  
PONE-D-16-20893 [pii]

Onofri, S., de Vera, J. P., Zucconi, L., Selbmann, L., Scalzi, G., Venkateswaran, K. J., Rabbow, E., de la Torre, R., and Horneck, G. (2015). [Survival of Antarctic Cryptoendolithic Fungi in Simulated Martian Conditions On Board the International Space Station](#). *Astrobiology* 15, 1052-1059. 10.1089/ast.2015.1324

Pacelli, C., Selbmann, L., Zucconi, L., De Vera, J. P., Rabbow, E., Horneck, G., de la Torre, R., and Onofri, S. (2016). [BIOMEX Experiment: Ultrastructural Alterations, Molecular Damage and Survival of the Fungus \*Cryomyces antarcticus\* after the Experiment Verification Tests](#). *Orig Life Evol Biosph.* 10.1007/s11084-016-9485-2

Tantikachornkiat, M., Sakakibara, S., Neuner, M., and Durall, D. M. (2016). [The use of propidium monoazide in conjunction with qPCR and Illumina sequencing to identify and quantify live yeasts and bacteria](#). *Int J Food Microbiol* 234, 53-59. 10.1016/j.ijfoodmicro.2016.06.031

Toledo Del Arbol, J., Perez Pulido, R., La Stora, A., Grande Burgos, M. J., Lucas, R., Ercolini, D., and Galvez, A. (2016). [Microbial diversity in pitted sweet cherries \(\*Prunus avium\* L.\) as affected by High-Hydrostatic Pressure treatment](#). *Food Res Int* 89, 790-796. S0963-9969(16)30435-5

Vendrame, M., Manzano, M., Comi, G., Bertrand, J., and Iacumin, L. (2014). [Use of propidium monoazide for the enumeration of viable \*Brettanomyces bruxellensis\* in wine and beer by quantitative PCR](#). *Food Microbiol* 42, 196-204. 10.1016/j.fm.2014.03.010

Venkateswaran, K., Vaishampayan, P., Cisneros, J., Pierson, D. L., Rogers, S. O., and Perry, J. (2014). [International Space Station environmental microbiome - microbial inventories of ISS filter debris](#). *Appl Microbiol Biotechnol* 98, 6453-6466. 10.1007/s00253-014-5650-6

Vesper, S., McKinstry, C., Hartmann, C., Neace, M., Yoder, S., and Vesper, A. (2008). [Quantifying fungal viability in air and water samples using quantitative PCR after treatment with propidium monoazide \(PMA\)](#). *J Microbiol Methods* 72, 180-184. 10.1016/j.mimet.2007.11.017

Wagner, A. O., Praeg, N., Reitschuler, C., and Illmer, P. (2015). [Effect of DNA extraction procedure, repeated extraction and ethidium monoazide \(EMA\)/propidium monoazide \(PMA\) treatment on overall DNA yield and impact on microbial fingerprints for bacteria, fungi and archaea in a reference soil](#). *Appl Soil Ecol* 93, 56-64. 10.1016/j.apsoil.2015.04.005

Weinmaier, T., Probst, A. J., La Duc, M. T., Ciobanu, D., Cheng, J. F., Ivanova, N., Rattei, T., and Vaishampayan, P. (2015). [A viability-linked metagenomic analysis of cleanroom environments: eukarya, prokaryotes, and viruses](#). *Microbiome* 3, 62. 10.1186/s40168-015-0129-y

## **Biofilms**

Alvarez, G., Gonzalez, M., Isabal, S., Blanc, V., and Leon, R. (2013). [Method to quantify live and dead cells in multi-species oral biofilm by real-time PCR with propidium monoazide](#). *AMB Express* 3, 1. 10.1186/2191-0855-3-1

de Almeida, J., Hoogenkamp, M., Felipe, W. T., Crielaard, W., and van der Waal, S. V. (2016). [Effectiveness of EDTA and Modified Salt Solution to Detach and Kill Cells from \*Enterococcus faecalis\* Biofilm](#). *J Endod* 42, 320-323. S0099-2399(15)01080-8

Fernandez Ramirez, M. D., Kostopoulos, I., Smid, E. J., Nierop Groot, M. N., and Abee, T. (2017). [Quantitative assessment of viable cells of \*Lactobacillus plantarum\* strains in single, dual and multi-strain biofilms](#). *Int J Food Microbiol* 244, 43-51. S0168-1605(16)30662-6

Fernandez, Y. M. M., Exterkate, R. A., Buijs, M. J., Crielaard, W., and Zaura, E. (2016). [Effect of mouthwashes on the composition and metabolic activity of oral biofilms grown in vitro](#). *Clin Oral Investig.* 10.1007/s00784-016-1876-2

Fernandez, Y. M. M., Exterkate, R. A. M., Buijs, M. J., Beertsen, W., van der Weijden, G. A., Zaura, E., and Crielaard, W. (2017). [A reproducible microcosm biofilm model of subgingival microbial communities](#). *J Periodontal Res* 52, 1021-1031. 10.1111/jre.12473

- Forbes, S., Cowley, N., Humphreys, G., Mistry, H., Amezcua, A., and McBain, A. J. (2017). [Formulation of Biocides Increases Antimicrobial Potency and Mitigates the Enrichment of Nonsusceptible Bacteria in Multispecies Biofilms](#). *Appl Environ Microbiol* 83. AEM.03054-16
- Gomez-Alvarez, V., Schrantz, K. A., Pressman, J. G., and Wahman, D. G. (2014). [Biofilm community dynamics in bench-scale annular reactors simulating arrestment of chloraminated drinking water nitrification](#). *Environ Sci Technol* 48, 5448-5457. 10.1021/es5005208
- Herrero, E. R., Slomka, V., Boon, N., Bernaerts, K., Hernandez-Sanabria, E., Quirynen, M., and Teughels, W. (2016). [Dysbiosis by neutralizing commensal mediated inhibition of pathobionts](#). *Sci Rep* 6, 38179. srep38179
- Kistler, J. O., Pesaro, M., and Wade, W. G. (2015). [Development and pyrosequencing analysis of an in-vitro oral biofilm model](#). *BMC Microbiol* 15, 24. 10.1186/s12866-015-0364-1
- Klein, M. I., Scott-Anne, K. M., Gregoire, S., Rosalen, P. L., and Koo, H. (2012). [Molecular approaches for viable bacterial population and transcriptional analyses in a rodent model of dental caries](#). *Mol Oral Microbiol* 27, 350-361. 10.1111/j.2041-1014.2012.00647.x
- Luo, Y., Bolt, H. L., Eggimann, G. A., McAuley, D. F., McMullan, R., Curran, T., Zhou, M., Jahoda, P. C., Cobb, S. L., and Lundy, F. T. (2016). [Peptoid Efficacy against Polymicrobial Biofilms Determined by Using Propidium Monoazide-Modified Quantitative PCR](#). *Chembiochem*. 10.1002/cbic.201600381
- Magajna, B., and Schraft, H. (2015). [Evaluation of Propidium Monoazide and Quantitative PCR To Quantify Viable \*Campylobacter jejuni\* Biofilm and Planktonic Cells in Log Phase and in a Viable but Nonculturable State](#). *J Food Prot* 78, 1303-1311. 10.4315/0362-028X.JFP-14-583
- Overney, A., Jacques-Andre-Coquin, J., Ng, P., Carpentier, B., Guillier, L., and Firmesse, O. (2017). [Impact of environmental factors on the culturability and viability of \*Listeria monocytogenes\* under conditions encountered in food processing plants](#). *Int J Food Microbiol* 244, 74-81. S0168-1605(16)30660-2
- Pan, Y., and Breidt, F., Jr. (2007). [Enumeration of viable \*Listeria monocytogenes\* cells by real-time PCR with propidium monoazide and ethidium monoazide in the presence of dead cells](#). *Appl Environ Microbiol* 73, 8028-8031. 10.1128/AEM.01198-07
- Pisz, J. M., Lawrence, J. R., Schafer, A. N., and Siciliano, S. D. (2007). [Differentiation of genes extracted from non-viable versus viable micro-organisms in environmental samples using ethidium monoazide bromide](#). *J Microbiol Methods* 71, 312-318. 10.1016/j.mimet.2007.09.015
- Sanchez, M. C., Marin, M. J., Figuero, E., Llama-Palacios, A., Leon, R., Blanc, V., Herrera, D., and Sanz, M. (2014). [Quantitative real-time PCR combined with propidium monoazide for the selective quantification of viable periodontal pathogens in an in vitro subgingival biofilm model](#). *J Periodontol Res* 49, 20-28. 10.1111/jre.12073
- Tavernier, S., and Coenye, T. (2015). [Quantification of \*Pseudomonas aeruginosa\* in multispecies biofilms using PMA-qPCR](#). *PeerJ* 3, e787. 10.7717/peerj.787
- Taylor, M. J., Bentham, R. H., and Ross, K. E. (2014). [Limitations of Using Propidium Monoazide with qPCR to Discriminate between Live and Dead \*Legionella\* in Biofilm Samples](#). *Microbiol Insights* 7, 15-24. 10.4137/MBI.S17723
- Vezzulli, L., Pezzati, E., Stauder, M., Stagnaro, L., Venier, P., and Pruzzo, C. (2014). [Aquatic ecology of the oyster pathogens \*Vibrio splendidus\* and \*Vibrio aestuarianus\*](#). *Environ Microbiol*. 10.1111/1462-2920.12484

## Eukaryotes

- Agullo-Barcelo, M., Moss, J. A., Green, J., Gillespie, S., Codony, F., Lucena, F., and Nocker, A. (2014). [Quantification of relative proportions of intact cells in microbiological samples using the example of \*Cryptosporidium parvum\* oocysts](#). *Lett Appl Microbiol* 58, 70-78. 10.1111/lam.12157
- Alonso, J. L., Amorós, I., and Guy, R. A. (2014). [Quantification of viable \*Giardia\* cysts and \*Cryptosporidium\* oocysts in wastewater using propidium monoazide quantitative real-time PCR](#). *Parasitol Res* 113, 2671-2678. 10.1007/s00436-014-3922-9
- Auld, R. R., Mykytczuk, N. C., Leduc, L. G., and Merritt, T. J. (2017). [Seasonal variation in an acid mine drainage microbial community](#). *Can J Microbiol* 63, 137-152. 10.1139/cjm-2016-0215
- Brescia, C. C., Griffin, S. M., Ware, M. W., Varughese, E. A., Egorov, A. I., and Villegas, E. N. (2009). [Cryptosporidium propidium monoazide-PCR, a molecular biology-based technique for genotyping of viable \*Cryptosporidium\* oocysts](#). *Appl Environ Microbiol* 75, 6856-6863. 10.1128/AEM.00540-09
- Burnet, J. B., Faraj, T., Cauchie, H. M., Joaquim-Justo, C., Servais, P., Prevost, M., and Dorner, S. M. (2017). [How does the cladoceran \*Daphnia pulex\* affect the fate of \*Escherichia coli\* in water?](#) *PLoS One* 12, e0171705.
- Fittipaldi, M., Pino Rodriguez, N. J., Adrados, B., Agusti, G., Penuela, G., Morato, J., and Codony, F. (2011). [Discrimination of viable \*Acanthamoeba castellanii\* trophozoites and cysts by propidium monoazide real-time polymerase chain reaction](#). *J Eukaryot Microbiol* 58, 359-364. 10.1111/j.1550-7408.2011.00557.x
- Habtewold, T., Groom, Z., Duchateau, L., and Christophides, G. K. (2015). [Detection of viable plasmodium ookinetes in the midguts of anopheles coluzzi using PMA-qrtPCR](#). *Parasit Vectors* 8, 455. 10.1186/s13071-015-1087-8
- Halfhide, C. P., Flanagan, B. F., Brearey, S. P., Hunt, J. A., Fonceca, A. M., McNamara, P. S., Howarth, D., Edwards, S., and Smyth, R. L. (2011). [Respiratory syncytial virus binds and undergoes transcription in neutrophils from the blood and airways of infants with severe bronchiolitis](#). *J Infect Dis* 204, 451-458. 10.1093/infdis/jir280
- Moreno-Mesonero, L., Moreno, Y., Alonso, J. L., and Ferrus, M. A. (2017). [Detection of viable \*Helicobacter pylori\* inside free-living amoebae in wastewater and drinking water samples from Eastern Spain](#). *Environ Microbiol* 19, 4103-4112. 10.1111/1462-2920.13856

## Viruses

- Bae, S., and Wuertz, S. (2012). [Survival of host-associated bacteroidales cells and their relationship with \*Enterococcus\* spp., \*Campylobacter jejuni\*, \*Salmonella enterica\* serovar Typhimurium, and adenovirus in freshwater microcosms as measured by propidium monoazide-quantitative PCR](#). *Appl Environ Microbiol* 78, 922-932. 10.1128/AEM.05157-11
- Bellehumeur, C., Boyle, B., Charette, S. J., Harel, J., L'Homme, Y., Masson, L., and Gagnon, C. A. (2015). [Propidium monoazide \(PMA\) and ethidium bromide monoazide \(EMA\) improve DNA array and high-throughput sequencing of porcine reproductive and respiratory syndrome virus identification](#). *J Virol Methods* 222, 182-191. 10.1016/j.jviromet.2015.06.014
- Bellehumeur, C., Nielsen, O., Measures, L., Harwood, L., Goldstein, T., Boyle, B., and Gagnon, C. A. (2016). [Herpesviruses including novel gammaherpesviruses are widespread among Phocid seal species in Canada](#). *J Wildl Dis* 52, 70-81. 10.7589/2015-01-020
- Corbeil, S., McColl, K. A., Williams, L. M., Slater, J., and Crane, M. S. J. (2017). [Innate resistance of New Zealand paua to abalone viral ganglioneuritis](#). *J Invertebr Pathol* 146, 31-35. S0022-2011(17)30081-2

- Coudray-Meunier, C.,Fraise, A.,Martin-Latil, S.,Guillier, L., and Perelle, S. (2013). [Discrimination of infectious hepatitis A virus and rotavirus by combining dyes and surfactants with RT-qPCR](#). BMC Microbiol 13, 216. 10.1186/1471-2180-13-216
- Elizaquivel, P.,Aznar, R., and Sanchez, G. (2013). [Recent developments in the use of viability dyes and quantitative PCR in the food microbiology field](#). J Appl Microbiol 116, 1-13. 10.1111/jam.12365
- Elsasser, D.,Ho, J.,Niessner, R.,Tiehm, A., and Seidel, M. (2018). [Heterogeneous asymmetric recombinase polymerase amplification \(haRPA\) for rapid hygiene control of large-volume water samples](#). Anal Biochem 546, 58-64. S0003-2697(18)30071-X
- Escudero-Abarca, B. I.,Rawsthorne, H.,Goulter, R. M.,Suh, S. H., and Jaykus, L. A. (2014). [Molecular methods used to estimate thermal inactivation of a prototype human norovirus: more heat resistant than previously believed?](#) Food Microbiol 41, 91-95. 10.1016/j.fm.2014.01.009
- Fittipaldi, M.,Rodriguez, N. J.,Codony, F.,Adrados, B.,Penuela, G. A., and Morato, J. (2010). [Discrimination of infectious bacteriophage T4 virus by propidium monoazide real-time PCR](#). J Virol Methods 168, 228-232. 10.1016/j.jviromet.2010.06.011
- Fraise, A.,Niveau, F.,Hennechart-Collette, C.,Coudray-Meunier, C.,Martin-Latil, S., and Perelle, S. (2018). [Discrimination of infectious and heat-treated norovirus by combining platinum compounds and real-time RT-PCR](#). Int J Food Microbiol 269, 64-74. S0168-1605(18)30023-0
- Garson, J. A.,Patel, P.,McDonald, C.,Ball, J.,Rosenberg, G.,Tettmar, K. I.,Brailsford, S. R.,Pitt, T., and Tedder, R. S. (2014). [Evaluation of an ethidium monoazide-enhanced 16S rDNA real-time polymerase chain reaction assay for bacterial screening of platelet concentrates and comparison with automated culture](#). Transfusion 54, 870-878. 10.1111/trf.12256
- Graiver, D. A.,Saunders, S. E.,Topliff, C. L.,Kelling, C. L., and Bartelt-Hunt, S. L. (2010). [Ethidium monoazide does not inhibit RT-PCR amplification of nonviable avian influenza RNA](#). J Virol Methods 164, 51-54. 10.1016/j.jviromet.2009.11.024
- Greening, G. (2013). [Foodborne viruses: a focus on challenges associated with detection methods](#). Microbiology Australia DOI: 10.1071/MA13022. 10.1071/MA13022
- Halfhide, C. P.,Flanagan, B. F.,Brearey, S. P.,Hunt, J. A.,Fonceca, A. M.,McNamara, P. S.,Howarth, D.,Edwards, S., and Smyth, R. L. (2011). [Respiratory syncytial virus binds and undergoes transcription in neutrophils from the blood and airways of infants with severe bronchiolitis](#). J Infect Dis 204, 451-458. jir280
- Karim, M. R.,Fout, G. S.,Johnson, C. H.,White, K. M., and Parshionkar, S. U. (2015). [Propidium monoazide reverse transcriptase PCR and RT-qPCR for detecting infectious enterovirus and norovirus](#). J Virol Methods 219, 51-61. S0166-0934(15)00042-7
- Kim, K.,Katayama, H.,Kitajima, M.,Tohya, Y., and Ohgaki, S. (2011). [Development of a real-time RT-PCR assay combined with ethidium monoazide treatment for RNA viruses and its application to detect viral RNA after heat exposure](#). Water Sci Technol 63, 502-507. 10.2166/wst.2011.249
- Kim, S. Y., and Ko, G. (2012). [Using propidium monoazide to distinguish between viable and nonviable bacteria, MS2 and murine norovirus](#). Lett Appl Microbiol 55, 182-188. 10.1111/j.1472-765X.2012.03276.x
- Lee, M.,Seo, D. J.,Seo, J.,Oh, H.,Jeon, S. B.,Ha, S. D.,Myoung, J.,Choi, I. S., and Choi, C. (2015). [Detection of viable murine norovirus using the plaque assay and propidium-monoazide-combined real-time reverse transcription-polymerase chain reaction](#). J Virol Methods. S0166-0934(15)00162-7

- Leifels, M., Hamza, I. A., Krieger, M., Wilhelm, M., Mackowiak, M., and Jurzik, L. (2016). [From Lab to Lake - Evaluation of Current Molecular Methods for the Detection of Infectious Enteric Viruses in Complex Water Matrices in an Urban Area](#). *PLoS One* 11, e0167105. 10.1371/journal.pone.0167105
- Leifels, M., Jurzik, L., Wilhelm, M., and Hamza, I. A. (2015). [Use of ethidium monoazide and propidium monoazide to determine viral infectivity upon inactivation by heat, UV- exposure and chlorine](#). *Int J Hyg Environ Health*. S1438-4639(15)00020-6
- Luo, Y., Bolt, H. L., Eggimann, G. A., McAuley, D. F., McMullan, R., Curran, T., Zhou, M., Jahoda, P. C., Cobb, S. L., and Lundy, F. T. (2016). [Peptoid Efficacy against Polymicrobial Biofilms Determined by Using Propidium Monoazide-Modified Quantitative PCR](#). *Chembiochem*. 10.1002/cbic.201600381
- McLellan, N. L., Lee, H., and Habash, M. B. (2016). [Evaluation of propidium monoazide and long-amplicon qPCR as an infectivity assay for coliphage](#). *J Virol Methods* 238, 48-55. S0166-0934(16)30086-6
- Moreno, L., Aznar, R., and Sanchez, G. (2015). [Application of viability PCR to discriminate the infectivity of hepatitis A virus in food samples](#). *Int J Food Microbiol* 201, 1-6. S0168-1605(15)00082-3
- Park, D., Shahbaz, H. M., Kim, S. H., Lee, M., Lee, W., Oh, J. W., Lee, D. U., and Park, J. (2016). [Inactivation efficiency and mechanism of UV-TiO<sub>2</sub> photocatalysis against murine norovirus using a solidified agar matrix](#). *Int J Food Microbiol* 238, 256-264. S0168-1605(16)30503-7
- Parshionikar, S., Laseke, I., and Fout, G. S. (2010). [Use of propidium monoazide in reverse transcriptase PCR to distinguish between infectious and noninfectious enteric viruses in water samples](#). *Appl Environ Microbiol* 76, 4318-4326. 10.1128/AEM.02800-09
- Prevost, B., Goulet, M., Lucas, F. S., Joyeux, M., Moulin, L., and Wurtzer, S. (2016). [Viral persistence in surface and drinking water: Suitability of PCR pre-treatment with intercalating dyes](#). *Water Res* 91, 68-76. 10.1016/j.watres.2015.12.049
- Quijada, N. M., Fongaro, G., Barardi, C. R., Hernandez, M., and Rodriguez-Lazaro, D. (2016). [Propidium Monoazide Integrated with qPCR Enables the Detection and Enumeration of Infectious Enteric RNA and DNA Viruses in Clam and Fermented Sausages](#). *Front Microbiol* 7, 2008. 10.3389/fmicb.2016.02008
- Randazzo, W., Lopez-Galvez, F., Allende, A., Aznar, R., and Sanchez, G. (2016). [Evaluation of viability PCR performance for assessing norovirus infectivity in fresh-cut vegetables and irrigation water](#). *Int J Food Microbiol* 229, 1-6. S0168-1605(16)30169-6
- Randazzo, W., Piqueras, J., Rodriguez-Diaz, J., Aznar, R., and Sanchez, G. (2018). [Improving efficiency of viability-qPCR for selective detection of infectious HAV in food and water samples](#). *J Appl Microbiol* 124, 958-964. 10.1111/jam.13519
- Sanchez, G., Elizaquivel, P., and Aznar, R. (2012). [Discrimination of infectious hepatitis A viruses by propidium monoazide real-time RT-PCR](#). *Food Environ Virol* 4, 21-25. 10.1007/s12560-011-9074-5
- Takahashi, H., Nakazawa, M., Ohshima, C., Sato, M., Tsuchiya, T., Takeuchi, A., Kunou, M., Kuda, T., and Kimura, B. (2015). [Heat-Denatured Lysozyme Inactivates Murine Norovirus as a Surrogate Human Norovirus](#). *Sci Rep* 5, 11819. 10.1038/srep11819
- Varma, M., Field, R., Stinson, M., Rukovets, B., Wymer, L., and Haugland, R. (2009). [Quantitative real-time PCR analysis of total and propidium monoazide-resistant fecal indicator bacteria in wastewater](#). *Water Res* 43, 4790-4801. 10.1016/j.watres.2009.05.031



Weinmaier, T., Probst, A. J., La Duc, M. T., Ciobanu, D., Cheng, J. F., Ivanova, N., Rattei, T., and Vaishampayan, P. (2015). [A viability-linked metagenomic analysis of cleanroom environments: eukarya, prokaryotes, and viruses](#). *Microbiome* 3, 62. 10.1186/s40168-015-0129-y

### Archaea

Auld, R. R., Mykytczuk, N. C., Leduc, L. G., and Merritt, T. J. (2017). [Seasonal variation in an acid mine drainage microbial community](#). *Can J Microbiol* 63, 137-152. 10.1139/cjm-2016-0215

Di Maiuta, N., Rufenacht, A., and Kuenzi, P. (2017). [Assessment of bacteria and archaea in metalworking fluids using massive parallel 16S rRNA gene tag sequencing](#). *Lett Appl Microbiol* 65, 266-273. 10.1111/lam.12782

Gagen, E. J., Huber, H., Meador, T., Hinrichs, K. U., and Thomm, M. (2013). [Novel cultivation-based approach to understanding the miscellaneous crenarchaeotic group \(MCG\) archaea from sedimentary ecosystems](#). *Appl Environ Microbiol* 79, 6400-6406. 10.1128/AEM.02153-13

Heise, J., Nega, M., Alawi, M., and Wagner, D. (2016). [Propidium monoazide treatment to distinguish between live and dead methanogens in pure cultures and environmental samples](#). *J Microbiol Methods* 121, 11-23. 10.1016/j.mimet.2015.12.002

Schirmack, J., Alawi, M., and Wagner, D. (2015). [Influence of Martian regolith analogs on the activity and growth of methanogenic archaea, with special regard to long-term desiccation](#). *Front Microbiol* 6, 210. 10.3389/fmicb.2015.00210

Wagner, A. O., Praeg, N., Reitschuler, C., and Illmer, P. (2015). [Effect of DNA extraction procedure, repeated extraction and ethidium monoazide \(EMA\)/propidium monoazide \(PMA\) treatment on overall DNA yield and impact on microbial fingerprints for bacteria, fungi and archaea in a reference soil](#). *Appl Soil Ecol* 93, 56-64. 10.1016/j.apsoil.2015.04.005

Weinmaier, T., Probst, A. J., La Duc, M. T., Ciobanu, D., Cheng, J. F., Ivanova, N., Rattei, T., and Vaishampayan, P. (2015). [A viability-linked metagenomic analysis of cleanroom environments: eukarya, prokaryotes, and viruses](#). *Microbiome* 3, 62. 10.1186/s40168-015-0129-y

Yergeau, E., Hogues, H., Whyte, L. G., and Greer, C. W. (2010). [The functional potential of high Arctic permafrost revealed by metagenomic sequencing, qPCR and microarray analyses](#). *ISME J* 4, 1206-1214. 10.1038/ismej.2010.41

### Food Science

Atia, A., Gomaa, A., Fernandez, B., Subirade, M., and Fliss, I. (2017). [Study and Understanding Behavior of Alginate-Inulin Synbiotics Beads for Protection and Delivery of Antimicrobial-Producing Probiotics in Colonic Simulated Conditions](#). *Probiotics Antimicrob Proteins*. 10.1007/s12602-017-9355-x

Caldwell, J. M., Juvonen, R., Brown, J., and Breidt, F. (2013). [Pectinatus sottacetoni sp. nov. isolated from commercial pickle spoilage tank](#). *Int J Syst Evol Microbiol* DOI: 10.1099/ijs.0.047886-0. 10.1099/ijs.0.047886-0

Cattani, F., Barth, V. C., Jr., Nasario, J. S., Ferreira, C. A., and Oliveira, S. D. (2016). [Detection and quantification of viable Bacillus cereus group species in milk by propidium monoazide quantitative real-time PCR](#). *J Dairy Sci*. 10.3168/jds.2015-10019

Cattani, F., Ferreira, C. A., and Oliveira, S. D. (2013). [The detection of viable vegetative cells of Bacillus sporothermodurans using propidium monoazide with semi-nested PCR](#). *Food Microbiol* 34, 196-201. 10.1016/j.fm.2012.12.007

- Crespo-Sempere, A., Estiarte, N., Marin, S., Sanchis, V., and Ramos, A. J. (2013). [Propidium monoazide combined with real-time quantitative PCR to quantify viable \*Alternaria\* spp. contamination in tomato products](#). *Int J Food Microbiol* 165, 214-220. 10.1016/j.ijfoodmicro.2013.05.017
- Desfosses-Foucault, E., Dussault-Lepage, V., Le Boucher, C., Savard, P., Lapointe, G., and Roy, D. (2012). [Assessment of Probiotic Viability during Cheddar Cheese Manufacture and Ripening Using Propidium Monoazide-PCR Quantification](#). *Front Microbiol* 3, 350. 10.3389/fmicb.2012.00350
- Duarte, A., Botteldoorn, N., Coucke, W., Denayer, S., Dierick, K., and Uyttendaele, M. (2015). [Effect of exposure to stress conditions on propidium monoazide \(PMA\)-qPCR based \*Campylobacter\* enumeration in broiler carcass rinses](#). *Food Microbiol* 48, 182-190. S0740-0020(14)00324-4
- Elizaquivel, P., Aznar, R., and Sanchez, G. (2013). [Recent developments in the use of viability dyes and quantitative PCR in the food microbiology field](#). *J Appl Microbiol* 116, 1-13. 10.1111/jam.12365
- Elizaquivel, P., Sanchez, G., Selma, M. V., and Aznar, R. (2012). [Application of propidium monoazide-qPCR to evaluate the ultrasonic inactivation of \*Escherichia coli\* O157:H7 in fresh-cut vegetable wash water](#). *Food Microbiol* 30, 316-320. 10.1016/j.fm.2011.10.008
- Erkus, O., de Jager, V. C., Geene, R. T., van Alen-Boerrigter, I., Hazelwood, L., van Hijum, S. A., Kleerebezem, M., and Smid, E. J. (2016). [Use of propidium monoazide for selective profiling of viable microbial cells during Gouda cheese ripening](#). *Int J Food Microbiol* 228, 1-9. S0168-1605(16)30142-8
- Feng, K., Hu, W., Jiang, A., Sarengaowa, Xu, Y., Zou, Y., Yang, L., and Wang, X. (2016). [A Dual Filtration-Based Multiplex PCR Method for Simultaneous Detection of Viable \*Escherichia coli\* O157:H7, \*Listeria monocytogenes\*, and \*Staphylococcus aureus\* on Fresh-Cut Cantaloupe](#). *PLoS One* 11, e0166874. 10.1371/journal.pone.0166874
- Forghani, F., Kim, J. B., and Oh, D. H. (2014). [Enterotoxigenic profiling of emetic toxin- and enterotoxin-producing \*Bacillus cereus\*, isolated from food, environmental, and clinical samples by multiplex PCR](#). *J Food Sci* 79, M2288-2293. 10.1111/1750-3841.12666
- Ganesan, B., Weimer, B. C., Pinzon, J., Dao Kong, N., Rompato, G., Brothersen, C., and McMahon, D. J. (2014). [Probiotic bacteria survive in Cheddar cheese and modify populations of other lactic acid bacteria](#). *J Appl Microbiol* 116, 1642-1656. 10.1111/jam.12482
- Grande Burgos, M. J., Lopez Aguayo, M. D. C., Perez Pulido, R., Galvez, A., and Lucas, R. (2017). [Analysis of the microbiota of refrigerated chopped parsley after treatments with a coating containing enterocin AS-48 or by high-hydrostatic pressure](#). *Food Res Int* 99, 91-97. S0963-9969(17)30203-X
- Greening, G. (2013). [Foodborne viruses: a focus on challenges associated with detection methods](#). *Microbiology Australia* DOI: 10.1071/MA13022. 10.1071/MA13022
- Josefsen, M. H., Lofstrom, C., Hansen, T. B., Christensen, L. S., Olsen, J. E., and Hoorfar, J. (2010). [Rapid quantification of viable \*Campylobacter\* bacteria on chicken carcasses, using real-time PCR and propidium monoazide treatment, as a tool for quantitative risk assessment](#). *Appl Environ Microbiol* 76, 5097-5104. 10.1128/AEM.00411-10
- Ju, W., Moyné, A. L., and Marco, M. L. (2016). [RNA-Based Detection Does not Accurately Enumerate Living \*Escherichia coli\* O157:H7 Cells on Plants](#). *Front Microbiol* 7, 223. 10.3389/fmicb.2016.00223
- Li, B., and Chen, J. Q. (2013). [Development of a sensitive and specific qPCR assay in conjunction with propidium monoazide for enhanced detection of live \*Salmonella\* spp. in food](#). *BMC Microbiol* 13, 273. 10.1186/1471-2180-13-273

- Liang, N., Dong, J., Luo, L., and Li, Y. (2011). [Detection of Viable Salmonella in Lettuce by Propidium Monoazide Real-Time PCR](#). *J Food Sci* 76, M234-237. 10.1111/j.1750-3841.2011.02123.x
- Liu, Y., and Mustapha, A. (2014). [Detection of viable Escherichia coli O157:H7 in ground beef by propidium monoazide real-time PCR](#). *Int J Food Microbiol* 170, 48-54. 10.1016/j.ijfoodmicro.2013.10.026
- Lo, R., Turner, M. S., Weeks, M., and Bansal, N. (2016). [Culture-independent bacterial community profiling of carbon dioxide treated raw milk](#). *Int J Food Microbiol* 233, 81-89. S0168-1605(16)30308-7
- Lo, R., Xue, T., Weeks, M., Turner, M. S., and Bansal, N. (2016). [Inhibition of bacterial growth in sweet cheese whey by carbon dioxide as determined by culture-independent community profiling](#). *Int J Food Microbiol* 217, 20-28. 10.1016/j.ijfoodmicro.2015.10.003
- Mace, S., Mamlouk, K., Chipchakova, S., Prevost, H., Joffraud, J. J., Dalgaard, P., Pilet, M. F., and Dousset, X. (2013). [Development of a Rapid Real-Time PCR Method as a Tool To Quantify Viable Photobacterium phosphoreum Bacteria in Salmon \(Salmo salar\) Steaks](#). *Appl Environ Microbiol* 79, 2612-2619. 10.1128/AEM.03677-12
- Medina, E., Perez-Diaz, I. M., Breidt, F., Hayes, J., Franco, W., Butz, N., and Azcarate-Peril, M. A. (2016). [Bacterial Ecology of Fermented Cucumber Rising pH Spoilage as Determined by Nonculture-Based Methods](#). *J Food Sci* 81, M121-129. 10.1111/1750-3841.13158
- Moyne, A. L., Harris, L. J., and Marco, M. L. (2013). [Assessments of total and viable Escherichia coli O157:H7 on field and laboratory grown lettuce](#). *PLoS One* 8, e70643. 10.1371/journal.pone.0070643
- Overney, A., Chassaing, D., Carpentier, B., Guillier, L., and Firmesse, O. (2016). [Development of synthetic media mimicking food soils to study the behaviour of Listeria monocytogenes on stainless steel surfaces](#). *Int J Food Microbiol* 238, 7-14. S0168-1605(16)30436-6
- Overney, A., Jacques-Andre-Coquin, J., Ng, P., Carpentier, B., Guillier, L., and Firmesse, O. (2017). [Impact of environmental factors on the culturability and viability of Listeria monocytogenes under conditions encountered in food processing plants](#). *Int J Food Microbiol* 244, 74-81. S0168-1605(16)30660-2 [pii] 10.1016/j.ijfoodmicro.2016.12.012
- Pacholewicz, E., Swart, A., Lipman, L. J., Wagenaar, J. A., Havelaar, A. H., and Duim, B. (2013). [Propidium monoazide does not fully inhibit the detection of dead Campylobacter on broiler chicken carcasses by qPCR](#). *J Microbiol Methods* 95, 32-38. S0167-7012(13)00178-4
- Quijada, N. M., Fongaro, G., Barardi, C. R., Hernandez, M., and Rodriguez-Lazaro, D. (2016). [Propidium Monoazide Integrated with qPCR Enables the Detection and Enumeration of Infectious Enteric RNA and DNA Viruses in Clam and Fermented Sausages](#). *Front Microbiol* 7, 2008. 10.3389/fmicb.2016.02008
- Randazzo, W., Lopez-Galvez, F., Allende, A., Aznar, R., and Sanchez, G. (2016). [Evaluation of viability PCR performance for assessing norovirus infectivity in fresh-cut vegetables and irrigation water](#). *Int J Food Microbiol* 229, 1-6. S0168-1605(16)30169-6
- Randazzo, W., Piqueras, J., Rodriguez-Diaz, J., Aznar, R., and Sanchez, G. (2018). [Improving efficiency of viability-qPCR for selective detection of infectious HAV in food and water samples](#). *J Appl Microbiol* 124, 958-964. 10.1111/jam.13519
- Rueckert, A., Ronimus, R. S., and Morgan, H. W. (2005). [Rapid differentiation and enumeration of the total, viable vegetative cell and spore content of thermophilic bacilli in milk powders with reference to Anoxybacillus flavithermus](#). *J Appl Microbiol* 99, 1246-1255. JAM2728

- Scariot, M. C., Venturelli, G. L., Prudencio, E. S., and Arisi, A. C. M. (2018). [Quantification of \*Lactobacillus paracasei\* viable cells in probiotic yoghurt by propidium monoazide combined with quantitative PCR](#). *Int J Food Microbiol* 264, 1-7. S0168-1605(17)30452-X
- Seliwiorstow, T., Duarte, A., Bare, J., Botteldoorn, N., Dierick, K., Uyttendaele, M., and De Zutter, L. (2015). [Comparison of sample types and analytical methods for the detection of highly campylobacter-colonized broiler flocks at different stages in the poultry meat production chain](#). *Foodborne Pathog Dis* 12, 399-405. 10.1089/fpd.2014.1894
- Soejima, T., Minami, J., and Iwatsuki, K. (2012). [Rapid propidium monoazide PCR assay for the exclusive detection of viable Enterobacteriaceae cells in pasteurized milk](#). *J Dairy Sci* 95, 3634-3642. 10.3168/jds.2012-5360
- Tian, Q., Feng, J. J., Hu, J., and Zhao, W. J. (2016). [Selective detection of viable seed-borne \*Acidovorax citrulli\* by real-time PCR with propidium monoazide](#). *Sci Rep* 6, 35457. srep35457
- Toledo Del Arbol, J., Perez Pulido, R., La Storia, A., Grande Burgos, M. J., Lucas, R., Ercolini, D., and Galvez, A. (2016). [Changes in microbial diversity of brined green asparagus upon treatment with high hydrostatic pressure](#). *Int J Food Microbiol* 216, 1-8. 10.1016/j.ijfoodmicro.2015.09.001
- Toledo Del Arbol, J., Perez Pulido, R., La Storia, A., Grande Burgos, M. J., Lucas, R., Ercolini, D., and Galvez, A. (2016). [Microbial diversity in pitted sweet cherries \(\*Prunus avium\* L.\) as affected by High-Hydrostatic Pressure treatment](#). *Food Res Int* 89, 790-796. S0963-9969(16)30435-5
- Udomsil, N., Chen, S., Rodtong, S., and Yongsawatdigul, J. (2016). [Quantification of viable bacterial starter cultures of \*Virgibacillus\* sp. and \*Tetragenococcus halophilus\* in fish sauce fermentation by real-time quantitative PCR](#). *Food Microbiol* 57, 54-62. S0740-0020(16)00005-8
- Vendrame, M., Iacumin, L., Manzano, M., and Comi, G. (2013). [Use of propidium monoazide for the enumeration of viable \*Oenococcus oeni\* in must and wine by quantitative PCR](#). *Food Microbiology* 35, 49-57. <http://dx.doi.org/10.1016/j.fm.2013.02.007>
- Villarreal, M. L., Padilha, M., Vieira, A. D., Franco, B. D., Martinez, R. C., and Saad, S. M. (2013). [Advantageous Direct Quantification of Viable Closely Related Probiotics in Petit-Suisse Cheeses under In Vitro Gastrointestinal Conditions by Propidium Monoazide - qPCR](#). *PLoS One* 8, e82102. 10.1371/journal.pone.0082102
- Wang, L., Li, P., Yang, Y., Xu, H., Aguilar, ZP, Xu, H., Yang, L, Xu, F, Lai, W, Xiong, Y, Wei, H (2013). [Development of an IMS-PMA-PCR assay with internal amplification control for rapid and sensitive detection of viable \*Escherichia coli\* O157:H7 in milk](#). *International Dairy Journal* DOI: 10.1016/j.idairyj.2013.07.006.
- Yang, Y., Xu, F., Xu, H., Aguilar, Z. P., Niu, R., Yuan, Y., Sun, J., You, X., Lai, W., Xiong, Y., et al. (2013). [Magnetic nano-beads based separation combined with propidium monoazide treatment and multiplex PCR assay for simultaneous detection of viable \*Salmonella\* Typhimurium, \*Escherichia coli\* O157:H7 and \*Listeria monocytogenes\* in food products](#). *Food Microbiol* 34, 418-424. 10.1016/j.fm.2013.01.004
- Youn, S. Y., Jeong, O. M., Choi, B. K., Jung, S. C., and Kang, M. S. (2016). [Application of loop-mediated isothermal amplification with propidium monoazide treatment to detect live \*Salmonella\* in chicken carcasses](#). *Poult Sci*. pew341
- Yu, S., Yan, L., Wu, X., Li, F., Wang, D., and Xu, H. (2017). [Multiplex PCR coupled with propidium monoazide for the detection of viable \*Cronobacter sakazakii\*, \*Bacillus cereus\*, and \*Salmonella\* spp. in milk and milk products](#). *J Dairy Sci* 100, 7874-7882. S0022-0302(17)30709-9

Zhang, Z.,Feng, L.,Xu, H.,Liu, C.,Shah, N. P., and Wei, H. (2016). [Detection of viable enterotoxin-producing \*Bacillus cereus\* and analysis of toxigenicity from ready-to-eat foods and infant formula milk powder by multiplex PCR](#). J Dairy Sci 99, 1047-1055. S0022-0302(15)00896-6

Zhang, Z.,Liu, H.,Lou, Y.,Xiao, L.,Liao, C.,Malakar, P. K.,Pan, Y., and Zhao, Y. (2015). [Quantifying viable \*Vibrio parahaemolyticus\* and \*Listeria monocytogenes\* simultaneously in raw shrimp](#). Appl Microbiol Biotechnol 99, 6451-6462. 10.1007/s00253-015-6715-x

Zhang, Z.,Liu, W.,Xu, H.,Aguilar, Z. P.,Shah, N. P., and Wei, H. (2015). [Propidium monoazide combined with real-time PCR for selective detection of viable \*Staphylococcus aureus\* in milk powder and meat products](#). J Dairy Sci. S0022-0302(15)00010-7

Zhou, B.,Chen, B.,Wu, X.,Li, F.,Yu, P.,Aguilar, Z. P.,Wei, H., and Xu, H. (2016). [A new application of a sodium deoxycholate-propidium monoazide-quantitative PCR assay for rapid and sensitive detection of viable \*Cronobacter sakazakii\* in powdered infant formula](#). J Dairy Sci 99, 9550-9559. S0022-0302(16)30721-4

Zhu, R.-G.,Li, T.-P.,Jia, Y.-F., and Song, L.-F. (2012). [Quantitative study of viable \*Vibrio parahaemolyticus\* cells in raw seafood using propidium monoazide in combination with quantitative PCR](#). Journal of Microbiological Methods DOI: 10.1016/j.mimet.2012.05.019. 10.1016/j.mimet.2012.05.019

## Probiotics

Atia, A.,Gomaa, A.,Fernandez, B.,Subirade, M., and Fliss, I. (2017). [Study and Understanding Behavior of Alginate-Inulin Synbiotics Beads for Protection and Delivery of Antimicrobial-Producing Probiotics in Colonic Simulated Conditions](#). Probiotics Antimicrob Proteins. 10.1007/s12602-017-9355-x

Desfosses-Foucault, E.,Dussault-Lepage, V.,Le Boucher, C.,Savard, P.,Lapointe, G., and Roy, D. (2012). [Assessment of Probiotic Viability during Cheddar Cheese Manufacture and Ripening Using Propidium Monoazide-PCR Quantification](#). Front Microbiol 3, 350. 10.3389/fmicb.2012.00350

Ganesan, B.,Weimer, B. C.,Pinzon, J.,Dao Kong, N.,Rompato, G.,Brothersen, C., and McMahon, D. J. (2014). [Probiotic bacteria survive in Cheddar cheese and modify populations of other lactic acid bacteria](#). J Appl Microbiol 116, 1642-1656. 10.1111/jam.12482

Khodaei, N.,Fernandez, B.,Fliss, I., and Karboune, S. (2016). [Digestibility and prebiotic properties of potato rhamnogalacturonan I polysaccharide and its galactose-rich oligosaccharides/oligomers](#). Carbohydr Polym 136, 1074-1084. 10.1016/j.carbpol.2015.09.106

Kiran, F.,Mokrani, M., and Osmanagaoglu, O. (2015). [Effect of Encapsulation on Viability of \*Pediococcus pentosaceus\* OZF During Its Passage Through the Gastrointestinal Tract Model](#). Curr Microbiol 71, 95-105. 10.1007/s00284-015-0832-8

Kramer, M.,Obermajer, N.,Bogovic Matijasic, B.,Rogelj, I., and Kmetec, V. (2009). [Quantification of live and dead probiotic bacteria in lyophilised product by real-time PCR and by flow cytometry](#). Appl Microbiol Biotechnol 84, 1137-1147. 10.1007/s00253-009-2068-7

Lai, C. H.,Wu, S. R.,Pang, J. C.,Ramireddy, L.,Chiang, Y. C.,Lin, C. K., and Tsen, H. Y. (2017). [Designing primers and evaluation of the efficiency of propidium monoazide - Quantitative polymerase chain reaction for counting the viable cells of \*Lactobacillus gasseri\* and \*Lactobacillus salivarius\*](#). J Food Drug Anal 25, 533-542. S1021-9498(16)30153-3

Oketic, K.,Bogovic Matijasic, B.,Obermajer, T.,Radulovic, Z.,Levic, S.,Mirkovic, N., and Nedovic, V. (2015). [Evaluation of propidium monoazide real-time PCR for enumeration of probiotic lactobacilli microencapsulated in calcium alginate beads](#). Benef Microbes, 1-9. C5406PK478061047

Scariot, M. C., Venturelli, G. L., Prudencio, E. S., and Arisi, A. C. M. (2018). [Quantification of \*Lactobacillus paracasei\* viable cells in probiotic yoghurt by propidium monoazide combined with quantitative PCR](#). *Int J Food Microbiol* 264, 1-7. S0168-1605(17)30452-X

Villarreal, M. L., Padilha, M., Vieira, A. D., Franco, B. D., Martinez, R. C., and Saad, S. M. (2013). [Advantageous Direct Quantification of Viable Closely Related Probiotics in Petit-Suisse Cheeses under In Vitro Gastrointestinal Conditions by Propidium Monoazide - qPCR](#). *PLoS One* 8, e82102. 10.1371/journal.pone.0082102

### **Environmental Testing**

Alonso, J. L., Amoros, I., and Guy, R. A. (2014). [Quantification of viable \*Giardia\* cysts and \*Cryptosporidium\* oocysts in wastewater using propidium monoazide quantitative real-time PCR](#). *Parasitol Res* 113, 2671-2678. 10.1007/s00436-014-3922-9

Bae, S., and Wuertz, S. (2009). [Rapid decay of host-specific fecal Bacteroidales cells in seawater as measured by quantitative PCR with propidium monoazide](#). *Water Res* 43, 4850-4859. 10.1016/j.watres.2009.06.053

Bae, S., and Wuertz, S. (2012). [Survival of host-associated bacteroidales cells and their relationship with \*Enterococcus\* spp., \*Campylobacter jejuni\*, \*Salmonella enterica\* serovar Typhimurium, and adenovirus in freshwater microcosms as measured by propidium monoazide-quantitative PCR](#). *Appl Environ Microbiol* 78, 922-932. 10.1128/AEM.05157-11

Bae, S., and Wuertz, S. (2014). [Decay of host-associated Bacteroidales cells and DNA in continuous-flow freshwater and seawater microcosms of identical experimental design and temperature as measured by PMA-qPCR and qPCR](#). *Water Res* 70C, 205-213. 10.1016/j.watres.2014.10.032

Chiao, T. H., Clancy, T. M., Pinto, A., Xi, C., and Raskin, L. (2014). [Differential resistance of drinking water bacterial populations to monochloramine disinfection](#). *Environ Sci Technol* 48, 4038-4047. 10.1021/es4055725

Desneux, J., Biscuit, A., Picard, S., and Pourcher, A. M. (2016). [Fate of Viable but Non-culturable \*Listeria monocytogenes\* in Pig Manure Microcosms](#). *Front Microbiol* 7, 245. 10.3389/fmicb.2016.00245

Desneux, J., Chemaly, M., and Pourcher, A. M. (2015). [Experimental design for the optimization of propidium monoazide treatment to quantify viable and non-viable bacteria in piggery effluents](#). *BMC Microbiol* 15, 164. 10.1186/s12866-015-0505-6

Ditommaso, S., Ricciardi, E., Giacomuzzi, M., Arauco Rivera, S. R., Ceccarelli, A., and Zotti, C. M. (2014). [Overestimation of the \*Legionella\* spp. load in environmental samples by quantitative real-time PCR: pretreatment with propidium monoazide as a tool for the assessment of an association between \*Legionella\* concentration and sanitary risk](#). *Diagn Microbiol Infect Dis* 80, 260-266. 10.1016/j.diagmicrobio.2014.09.010

Ditommaso, S., Ricciardi, E., Giacomuzzi, M., Arauco Rivera, S. R., and Zotti, C. M. (2014). [\*Legionella\* in water samples: How can you interpret the results obtained by quantitative PCR?](#) *Mol Cell Probes*. 10.1016/j.mcp.2014.09.002

Dong, S., Hong, P. Y., and Nguyen, T. H. (2014). [Persistence of \*Bacteroides ovatus\* under simulated sunlight irradiation](#). *BMC Microbiol* 14, 178. 10.1186/1471-2180-14-178

Eichmiller, J. J., Borchert, A. J., Sadowsky, M. J., and Hicks, R. E. (2014). [Decay of genetic markers for fecal bacterial indicators and pathogens in sand from Lake Superior](#). *Water Res* 59, 99-111. 10.1016/j.watres.2014.04.005

Elsasser, D., Ho, J., Niessner, R., Tiehm, A., and Seidel, M. (2018). [Heterogeneous asymmetric recombinase polymerase amplification \(haRPA\) for rapid hygiene control of large-volume water samples](#). *Anal Biochem* 546, 58-64. 10.1016/j.ab.2018.01.032

- Fittipaldi, M., Codony, F., Adrados, B., Camper, A. K., and Morato, J. (2010). [Viable real-time PCR in environmental samples: can all data be interpreted directly?](#) *Microb Ecol* 61, 7-12. 10.1007/s00248-010-9719-1
- Forghani, F., Kim, J. B., and Oh, D. H. (2014). [Enterotoxigenic profiling of emetic toxin- and enterotoxin-producing \*Bacillus cereus\*, Isolated from food, environmental, and clinical samples by multiplex PCR.](#) *J Food Sci* 79, M2288-2293. 10.1111/1750-3841.12666
- Fujimoto, M., Moyerbrailean, G. A., Noman, S., Gizicki, J. P., Ram, M. L., Green, P. A., and Ram, J. L. (2014). [Application of ion torrent sequencing to the assessment of the effect of alkali ballast water treatment on microbial community diversity.](#) *PLoS One* 9, e107534. 10.1371/journal.pone.0107534
- Gensberger, E. T., Polt, M., Konrad-Koszler, M., Kinner, P., Sessitsch, A., and Kostic, T. (2014). [Evaluation of quantitative PCR combined with PMA treatment for molecular assessment of microbial water quality.](#) *Water Res* 67, 367-376. 10.1016/j.watres.2014.09.022
- Heise, J., Nega, M., Alawi, M., and Wagner, D. (2016). [Propidium monoazide treatment to distinguish between live and dead methanogens in pure cultures and environmental samples.](#) *J Microbiol Methods* 121, 11-23. 10.1016/j.mimet.2015.12.002
- Kibbee, R. J., and Ormeci, B. (2016). [Development of a sensitive and false-positive free PMA-qPCR viability assay to quantify VBNC \*Escherichia coli\* and evaluate disinfection performance in wastewater effluent.](#) *J Microbiol Methods* 132, 139-147. 10.1016/j.mimet.2016.12.004
- Kim, M., and Wuertz, S. (2015). [Survival and persistence of host-associated Bacteroidales cells and DNA in comparison with \*Escherichia coli\* and \*Enterococcus\* in freshwater sediments as quantified by PMA-qPCR and qPCR.](#) *Water Res* 87, 182-192. 10.1016/j.watres.2015.09.014
- Leifels, M., Hamza, I. A., Krieger, M., Wilhelm, M., Mackowiak, M., and Jurzik, L. (2016). [From Lab to Lake - Evaluation of Current Molecular Methods for the Detection of Infectious Enteric Viruses in Complex Water Matrices in an Urban Area.](#) *PLoS One* 11, e0167105. 10.1371/journal.pone.0167105
- Li, R., Tun, H. M., Jahan, M., Zhang, Z., Kumar, A., Fernando, D., Farenhorst, A., and Khafipour, E. (2017). [Comparison of DNA-, PMA-, and RNA-based 16S rRNA Illumina sequencing for detection of live bacteria in water.](#) *Sci Rep* 7, 5752. 10.1038/s41598-017-02516-3
- Lu, C., Li, S., Gong, S., Yuan, S., and Yu, X. (2015). [Mixing regime as a key factor to determine DON formation in drinking water biological treatment.](#) *Chemosphere*. 10.1016/j.chemosphere.2014.12.059
- Magiopoulos, I., McQuillan, J. S., Burd, C. L., Mowlem, M., and Tsaloglou, M. N. (2016). [A multi-parametric assessment of decontamination protocols for the subglacial Lake Ellsworth probe.](#) *J Microbiol Methods*. 10.1016/j.mimet.2016.02.012
- Mayer, T., Blachowicz, A., Probst, A. J., Vaishampayan, P., Checinska, A., Swarmer, T., de Leon, P., and Venkateswaran, K. (2016). [Microbial succession in an inflated lunar/Mars analog habitat during a 30-day human occupation.](#) *Microbiome* 4, 22. 10.1186/s40168-016-0167-0
- Moreno-Mesonero, L., Moreno, Y., Alonso, J. L., and Ferrus, M. A. (2017). [Detection of viable \*Helicobacter pylori\* inside free-living amoebae in wastewater and drinking water samples from Eastern Spain.](#) *Environ Microbiol* 19, 4103-4112. 10.1111/1462-2920.13856
- Pang, Y. C., Xi, J. Y., Xu, Y., Huo, Z. Y., and Hu, H. Y. (2016). [Shifts of live bacterial community in secondary effluent by chlorine disinfection revealed by Miseq high-throughput sequencing combined with propidium monoazide treatment.](#) *Appl Microbiol Biotechnol* 100, 6435-6446. 10.1007/s00253-016-7452-5

- Parshionikar, S., Laseke, I., and Fout, G. S. (2010). [Use of propidium monoazide in reverse transcriptase PCR to distinguish between infectious and noninfectious enteric viruses in water samples](#). *Appl Environ Microbiol* 76, 4318-4326. 10.1128/AEM.02800-09
- Prevost, B., Goulet, M., Lucas, F. S., Joyeux, M., Moulin, L., and Wurtzer, S. (2016). [Viral persistence in surface and drinking water: Suitability of PCR pre-treatment with intercalating dyes](#). *Water Res* 91, 68-76. 10.1016/j.watres.2015.12.049
- Salam, K. W., El-Fadel, M., Barbour, E. K., and Saikaly, P. E. (2014). [A propidium monoazide-quantitative PCR method for the detection and quantification of viable \*Enterococcus faecalis\* in large-volume samples of marine waters](#). *Appl Microbiol Biotechnol*. 10.1007/s00253-014-6023-x
- Santiago, P., Moreno, Y., and Ferrus, M. A. (2015). [Identification of Viable \*Helicobacter pylori\* in Drinking Water Supplies by Cultural and Molecular Techniques](#). *Helicobacter*. 10.1111/hel.12205
- Scaturro, M., Fontana, S., Dell'eva, I., Helfer, F., Marchio, M., Stefanetti, M. V., Cavallaro, M., Miglietta, M., Montagna, M. T., De Giglio, O., et al. (2016). [A multicenter study of viable PCR using propidium monoazide to detect \*Legionella\* in water samples](#). *Diagn Microbiol Infect Dis* 85, 283-288. 10.1016/j.diagmicrobio.2016.04.009
- Singh, G., Vajpayee, P., Bhatti, S., Ronnie, N., Shah, N., McClure, P., and Shanker, R. (2013). [Determination of viable \*Salmonellae\* from potable and source water through PMA assisted qPCR](#). *Ecotoxicol Environ Saf*. 10.1016/j.ecoenv.2013.02.017
- Soto-Munoz, L., Teixido, N., Usall, J., Vinas, I., Crespo-Sempere, A., and Torres, R. (2014). [Development of PMA real-time PCR method to quantify viable cells of \*Pantoea agglomerans\* CPA-2, an antagonist to control the major postharvest diseases on oranges](#). *Int J Food Microbiol* 180, 49-55. 10.1016/j.ijfoodmicro.2014.04.011
- Tian, Q., Feng, J. J., Hu, J., and Zhao, W. J. (2016). [Selective detection of viable seed-borne \*Acidovorax citrulli\* by real-time PCR with propidium monoazide](#). *Sci Rep* 6, 35457. 10.1038/srep35457
- Truchado, P., Hernandez, N., Gil, M. I., Ivanek, R., and Allende, A. (2018). [Correlation between \*E. coli\* levels and the presence of foodborne pathogens in surface irrigation water: Establishment of a sampling program](#). *Water Res* 128, 226-233. 10.1016/j.watres.2017.10.041
- van Frankenhuyzen, J. K., Trevors, J. T., Flemming, C. A., Lee, H., and Habash, M. B. (2013). [Optimization, validation, and application of a real-time PCR protocol for quantification of viable bacterial cells in municipal sewage sludge and biosolids using reporter genes and \*Escherichia coli\*](#). *J Ind Microbiol Biotechnol* 40, 1251-1261. 10.1007/s10295-013-1319-x
- Varma, M., Field, R., Stinson, M., Rukovets, B., Wymer, L., and Haugland, R. (2009). [Quantitative real-time PCR analysis of total and propidium monoazide-resistant fecal indicator bacteria in wastewater](#). *Water Res* 43, 4790-4801. 10.1016/j.watres.2009.05.031
- Wicaksono, W. A., Jones, E. E., Monk, J., and Ridgway, H. J. (2016). [The Bacterial Signature of \*Leptospermum scoparium\* \(Manuka\) Reveals Core and Accessory Communities with Bioactive Properties](#). *PLoS One* 11, e0163717. 10.1371/journal.pone.0163717
- Yang, Y., Cheng, D., Li, Y., Yu, L., Gin, K. Y., Chen, J. P., and Reinhard, M. (2017). [Effects of monochloramine and hydrogen peroxide on the bacterial community shifts in biologically treated wastewater](#). *Chemosphere* 189, 399-406. 10.1016/j.chemosphere.2017.09.087



## Filtered Samples

Chiao, T. H., Clancy, T. M., Pinto, A., Xi, C., and Raskin, L. (2014). [Differential resistance of drinking water bacterial populations to monochloramine disinfection](#). *Environ Sci Technol* 48, 4038-4047. 10.1021/es4055725

Hellein, K. N., Kennedy, E. M., Harwood, V. J., Gordon, K. V., Wang, S. Y., and Lepo, J. E. (2012). [A filter-based propidium monoazide technique to distinguish live from membrane-compromised microorganisms using quantitative PCR](#). *J Microbiol Methods*. 10.1016/j.mimet.2012.01.015

Salam, K. W., El-Fadel, M., Barbour, E. K., and Saikaly, P. E. (2014). [A propidium monoazide-quantitative PCR method for the detection and quantification of viable \*Enterococcus faecalis\* in large-volume samples of marine waters](#). *Appl Microbiol Biotechnol*. 10.1007/s00253-014-6023-x

Slimani, S., Robyns, A., Jarraud, S., Molmeret, M., Dusserre, E., Mazure, C., Facon, J. P., Lina, G., Etienne, J., and Ginevra, C. (2012). [Evaluation of propidium monoazide \(PMA\) treatment directly on membrane filter for the enumeration of viable but non cultivable \*Legionella\* by qPCR](#). *J Microbiol Methods* 88, 319-321. 10.1016/j.mimet.2011.12.010

Tseng, C. C., Hsiao, P. K., Chang, K. C., Cheng, C. C., Yiin, L. M., and Hsieh, C. J. (2014). [Detection of Viable Antibiotic-Resistant/Sensitive \*Acinetobacter baumannii\* in Indoor Air by Propidium Monoazide Quantitative PCR](#). *Indoor Air*. 10.1111/ina.12165

Venkateswaran, K., Vaishampayan, P., Cisneros, J., Pierson, D. L., Rogers, S. O., and Perry, J. (2014). [International Space Station environmental microbiome - microbial inventories of ISS filter debris](#). *Appl Microbiol Biotechnol* 98, 6453-6466. 10.1007/s00253-014-5650-6

Vesper, S., McKinstry, C., Hartmann, C., Neace, M., Yoder, S., and Vesper, A. (2008). [Quantifying fungal viability in air and water samples using quantitative PCR after treatment with propidium monoazide \(PMA\)](#). *J Microbiol Methods* 72, 180-184. 10.1016/j.mimet.2007.11.017

## Sequencing

Auld, R. R., Mykytczuk, N. C., Leduc, L. G., and Merritt, T. J. (2017). [Seasonal variation in an acid mine drainage microbial community](#). *Can J Microbiol* 63, 137-152. 10.1139/cjm-2016-0215

Be, N. A., Avila-Herrera, A., Allen, J. E., Singh, N., Checinska Sielaff, A., Jaing, C., and Venkateswaran, K. (2017). [Whole metagenome profiles of particulates collected from the International Space Station](#). *Microbiome* 5, 81. 10.1186/s40168-017-0292-4

Bellehumeur, C., Boyle, B., Charette, S. J., Harel, J., L'Homme, Y., Masson, L., and Gagnon, C. A. (2015). [Propidium monoazide \(PMA\) and ethidium bromide monoazide \(EMA\) improve DNA array and high-throughput sequencing of porcine reproductive and respiratory syndrome virus identification](#). *J Virol Methods* 222, 182-191. 10.1016/j.jviromet.2015.06.014

Blachowicz, A., Mayer, T., Bashir, M., Pieber, T. R., De Leon, P., and Venkateswaran, K. (2017). [Human presence impacts fungal diversity of inflated lunar/Mars analog habitat](#). *Microbiome* 5, 62. 10.1186/s40168-017-0280-8

Boutin, S., Depner, M., Stahl, M., Graeber, S. Y., Dittrich, S. A., Legatzki, A., von Mutius, E., Mall, M., and Dalpke, A. H. (2017). [Comparison of Oropharyngeal Microbiota from Children with Asthma and Cystic Fibrosis](#). *Mediators Inflamm* 2017, 5047403. 10.1155/2017/5047403

Boutin, S., Graeber, S. Y., Weitnauer, M., Panitz, J., Stahl, M., Clauszntzer, D., Kaderali, L., Einarsson, G., Tunney, M. M., Elborn, J. S., et al. (2015). [Comparison of microbiomes from different niches of upper and lower airways in children and adolescents with cystic fibrosis](#). *PLoS One* 10, e0116029. 10.1371/journal.pone.0116029

- Checinska, A., Probst, A. J., Vaishampayan, P., White, J. R., Kumar, D., Stepanov, V. G., Fox, G. E., Nilsson, H. R., Pierson, D. L., Perry, J., et al. (2015). [Microbiomes of the dust particles collected from the International Space Station and Spacecraft Assembly Facilities](#). *Microbiome* 3, 50. 10.1186/s40168-015-0116-3
- Chiao, T. H., Clancy, T. M., Pinto, A., Xi, C., and Raskin, L. (2014). [Differential resistance of drinking water bacterial populations to monochloramine disinfection](#). *Environ Sci Technol* 48, 4038-4047. 10.1021/es4055725
- Cuthbertson, L., Rogers, G. B., Walker, A. W., Oliver, A., Green, L. E., Daniels, T. W., Carroll, M. P., Parkhill, J., Bruce, K. D., and van der Gast, C. J. (2015). [Respiratory microbiota resistance and resilience to pulmonary exacerbation and subsequent antimicrobial intervention](#). *ISME J.* 10.1038/ismej.2015.198
- Cuthbertson, L., Rogers, G. B., Walker, A. W., Oliver, A., Hafiz, T., Hoffman, L. R., Carroll, M. P., Parkhill, J., Bruce, K. D., and van der Gast, C. J. (2014). [Time between collection and storage significantly influences bacterial sequence composition in sputum samples from cystic fibrosis respiratory infections](#). *J Clin Microbiol* 52, 3011-3016. 10.1128/JCM.00764-14
- Cuthbertson, L., Rogers, G. B., Walker, A. W., Oliver, A., Hoffman, L. R., Carroll, M. P., Parkhill, J., Bruce, K. D., and van der Gast, C. J. (2014). [Implications of multiple freeze-thawing on respiratory samples for culture-independent analyses](#). *J Cyst Fibros.* S1569-1993(14)00231-8
- Di Maiuta, N., Rufenacht, A., and Kuenzi, P. (2017). [Assessment of bacteria and archaea in metalworking fluids using massive parallel 16S rRNA gene tag sequencing](#). *Lett Appl Microbiol* 65, 266-273. 10.1111/lam.12782
- Eramo, A., Medina, W. M., and Fahrenfeld, N. L. (2017). [Peracetic acid disinfection kinetics for combined sewer overflows: indicator organisms, antibiotic resistance genes, and microbial community](#). *Environ Sci (Camb)* 3, 1061-1072. 10.1039/C7EW00184C
- Erkus, O., de Jager, V. C., Geene, R. T., van Alen-Boerrigter, I., Hazelwood, L., van Hijum, S. A., Kleerebezem, M., and Smid, E. J. (2016). [Use of propidium monoazide for selective profiling of viable microbial cells during Gouda cheese ripening](#). *Int J Food Microbiol* 228, 1-9. S0168-1605(16)30142-8
- Exterkate, R. A., Zaura, E., Brandt, B. W., Buijs, M. J., Koopman, J. E., Crielaard, W., and Ten Cate, J. M. (2014). [The effect of propidium monoazide treatment on the measured bacterial composition of clinical samples after the use of a mouthwash](#). *Clin Oral Investig.* 10.1007/s00784-014-1297-z
- Fernandez, Y. M. M., Exterkate, R. A., Buijs, M. J., Crielaard, W., and Zaura, E. (2016). [Effect of mouthwashes on the composition and metabolic activity of oral biofilms grown in vitro](#). *Clin Oral Investig.* 10.1007/s00784-016-1876-2
- Fujimoto, M., Moyerbrailean, G. A., Noman, S., Gizicki, J. P., Ram, M. L., Green, P. A., and Ram, J. L. (2014). [Application of ion torrent sequencing to the assessment of the effect of alkali ballast water treatment on microbial community diversity](#). *PLoS One* 9, e107534. 10.1371/journal.pone.0107534
- Gomez-Alvarez, V., Schrantz, K. A., Pressman, J. G., and Wahman, D. G. (2014). [Biofilm community dynamics in bench-scale annular reactors simulating arrestment of chloraminated drinking water nitrification](#). *Environ Sci Technol* 48, 5448-5457. 10.1021/es5005208
- Guo, F., and Zhang, T. (2014). [Detecting the Nonviable and Heat-Tolerant Bacteria in Activated Sludge by Minimizing DNA from Dead Cells](#). *Microb Ecol.* 10.1007/s00248-014-0389-2
- Kistler, J. O., Pesaro, M., and Wade, W. G. (2015). [Development and pyrosequencing analysis of an in-vitro oral biofilm model](#). *BMC Microbiol* 15, 24. 10.1186/s12866-015-0364-1

- Li, R., Tun, H. M., Jahan, M., Zhang, Z., Kumar, A., Fernando, D., Farenhorst, A., and Khafipour, E. (2017). [Comparison of DNA-, PMA-, and RNA-based 16S rRNA Illumina sequencing for detection of live bacteria in water](#). *Sci Rep* 7, 5752. 10.1038/s41598-017-02516-3
- Li, Y. F., Chen, P. H., and Yu, Z. (2014). [Spatial and temporal variations of microbial community in a mixed plug-flow loop reactor fed with dairy manure](#). *Microb Biotechnol* 7, 332-346. 10.1111/1751-7915.12125
- Lo, R., Turner, M. S., Weeks, M., and Bansal, N. (2016). [Culture-independent bacterial community profiling of carbon dioxide treated raw milk](#). *Int J Food Microbiol* 233, 81-89. S0168-1605(16)30308-7
- Mahnert, A., Vaishampayan, P., Probst, A. J., Auerbach, A., Moissl-Eichinger, C., Venkateswaran, K., and Berg, G. (2015). [Cleanroom Maintenance Significantly Reduces Abundance but Not Diversity of Indoor Microbiomes](#). *PLoS One* 10, e0134848. 10.1371/journal.pone.0134848
- Mayer, T., Blachowicz, A., Probst, A. J., Vaishampayan, P., Checinska, A., Swarmer, T., de Leon, P., and Venkateswaran, K. (2016). [Microbial succession in an inflated lunar/Mars analog habitat during a 30-day human occupation](#). *Microbiome* 4, 22. 10.1186/s40168-016-0167-0
- Moissl-Eichinger, C., Auerbach, A. K., Probst, A. J., Mahnert, A., Tom, L., Piceno, Y., Andersen, G. L., Venkateswaran, K., Rettberg, P., Barczyk, S., et al. (2015). [Quo vadis? Microbial profiling revealed strong effects of cleanroom maintenance and routes of contamination in indoor environments](#). *Sci Rep* 5, 9156. srep09156
- Moreno-Mesonero, L., Moreno, Y., Alonso, J. L., and Ferrus, M. A. (2017). [Detection of viable \*Helicobacter pylori\* inside free-living amoebae in wastewater and drinking water samples from Eastern Spain](#). *Environ Microbiol* 19, 4103-4112. 10.1111/1462-2920.13856
- Nguyen, L. D., Deschaght, P., Merlin, S., Loywick, A., Audebert, C., Van Daele, S., Viscogliosi, E., Vaneechoutte, M., and Delhaes, L. (2016). [Effects of Propidium Monoazide \(PMA\) Treatment on Mycobiome and Bacteriome Analysis of Cystic Fibrosis Airways during Exacerbation](#). *PLoS One* 11, e0168860. 10.1371/journal.pone.0168860
- Nocker, A., Richter-Heitmann, T., Montijn, R., Schuren, F., and Kort, R. (2010). [Discrimination between live and dead cells in bacterial communities from environmental water samples analyzed by 454 pyrosequencing](#). *Int Microbiol* 13, 59-65
- Pang, Y. C., Xi, J. Y., Xu, Y., Huo, Z. Y., and Hu, H. Y. (2016). [Shifts of live bacterial community in secondary effluent by chlorine disinfection revealed by Miseq high-throughput sequencing combined with propidium monoazide treatment](#). *Appl Microbiol Biotechnol* 100, 6435-6446. 10.1007/s00253-016-7452-5
- Rogers, G. B., Cuthbertson, L., Hoffman, L. R., Wing, P. A., Pope, C., Hooftman, D. A., Lilley, A. K., Oliver, A., Carroll, M. P., Bruce, K. D., et al. (2012). [Reducing bias in bacterial community analysis of lower respiratory infections](#). *ISME J* DOI: 10.1038/ismej.2012.145. 10.1038/ismej.2012.145
- Rogers, G. B., van der Gast, C. J., Bruce, K. D., Marsh, P., Collins, J. E., Sutton, J., and Wright, M. (2013). [Ascitic microbiota composition is correlated with clinical severity in cirrhosis with portal hypertension](#). *PLoS One* 8, e74884. 10.1371/journal.pone.0074884
- Tantikachornkiat, M., Sakakibara, S., Neuner, M., and Durall, D. M. (2016). [The use of propidium monoazide in conjunction with qPCR and Illumina sequencing to identify and quantify live yeasts and bacteria](#). *Int J Food Microbiol* 234, 53-59. 10.1016/j.ijfoodmicro.2016.06.031
- Thompson, H., Rybalka, A., Moazzez, R., Dewhirst, F. E., and Wade, W. G. (2015). [In vitro culture of previously uncultured oral bacterial phylotypes](#). *Appl Environ Microbiol* 81, 8307-8314. 10.1128/AEM.02156-15

Toledo Del Arbol, J., Perez Pulido, R., La Stora, A., Grande Burgos, M. J., Lucas, R., Ercolini, D., and Galvez, A. (2016). [Changes in microbial diversity of brined green asparagus upon treatment with high hydrostatic pressure](#). *Int J Food Microbiol* 216, 1-8. 10.1016/j.ijfoodmicro.2015.09.001

Venkateswaran, K., Vaishampayan, P., Cisneros, J., Pierson, D. L., Rogers, S. O., and Perry, J. (2014). [International Space Station environmental microbiome - microbial inventories of ISS filter debris](#). *Appl Microbiol Biotechnol* 98, 6453-6466. 10.1007/s00253-014-5650-6

Weinmaier, T., Probst, A. J., La Duc, M. T., Ciobanu, D., Cheng, J. F., Ivanova, N., Rattei, T., and Vaishampayan, P. (2015). [A viability-linked metagenomic analysis of cleanroom environments: eukarya, prokaryotes, and viruses](#). *Microbiome* 3, 62. 10.1186/s40168-015-0129-y

Yang, Y., Cheng, D., Li, Y., Yu, L., Gin, K. Y., Chen, J. P., and Reinhard, M. (2017). [Effects of monochloramine and hydrogen peroxide on the bacterial community shifts in biologically treated wastewater](#). *Chemosphere* 189, 399-406. S0045-6535(17)31500-X

Yergeau, E., Hogues, H., Whyte, L. G., and Greer, C. W. (2010). [The functional potential of high Arctic permafrost revealed by metagenomic sequencing, qPCR and microarray analyses](#). *ISME J* 4, 1206-1214. 10.1038/ismej.2010.41

### **Loop-Mediated Isothermal Amplification (LAMP)**

Chen, S., Wang, F., Beaulieu, J. C., Stein, R. E., and Ge, B. (2011). [Rapid detection of viable salmonellae in produce by coupling propidium monoazide with loop-mediated isothermal amplification](#). *Appl Environ Microbiol* 77, 4008-4016. 10.1128/AEM.00354-11

Li, Y., Yang, L., Fu, J., Yan, M., Chen, D., and Zhang, L. (2017). [The novel loop-mediated isothermal amplification based confirmation methodology on the bacteria in Viable but Non-Culturable \(VBNC\) state](#). *Microb Pathog* 111, 280-284. S0882-4010(17)30999-3

Wan, C., Yang, Y., Xu, H., Aguilar, Z. P., Liu, C., Lai, W., Xiong, Y., Xu, F., and Wei, H. (2012). [Development of a propidium monoazide treatment combined with loop-mediated isothermal amplification \(PMA-LAMP\) assay for rapid detection of viable \*Listeria monocytogenes\*](#). *International Journal of Food Science & Technology* 47, 2460-2467. 10.1111/j.1365-2621.2012.03123.x

Yan, M., Xu, L., Jiang, H., Zhou, Z., Zhou, S., and Zhang, L. (2017). [PMA-LAMP for rapid detection of \*Escherichia coli\* and shiga toxins from viable but non-culturable state](#). *Microb Pathog* 105, 245-250. S0882-4010(16)30933-0 [pii] 10.1016/j.micpath.2017.02.001

Youn, S. Y., Jeong, O. M., Choi, B. K., Jung, S. C., and Kang, M. S. (2016). [Application of loop-mediated isothermal amplification with propidium monoazide treatment to detect live \*Salmonella\* in chicken carcasses](#). *Poult Sci*. pew341

Zhao, X., Wang, J., Forghani, F., Park, J. H., Park, M. S., Seo, K. H., and Oh, D. H. (2013). [Rapid Detection of Viable \*Escherichia coli\* O157 by Coupling Propidium Monoazide with Loop-Mediated Isothermal Amplification](#). *J Microbiol Biotechnol* 23, 1708-1716. 10.4014/jmb.1306.06003

Zhong, H., Zhong, Y., Deng, Q., Zhou, Z., Guan, X., Yan, M., Hu, T., and Luo, M. (2017). [Virulence of thermolabile haemolysin, gastroenteritis related pathogenicity toxin and toxin of the pathogens \*Vibrio Parahaemolyticus\* in Viable but Non-Culturable \(VBNC\) state](#). *Microb Pathog* 111, 352-356. S0882-4010(17)31068-9