
Selected EvaGreen® References by Application

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

Contents:

The use of EvaGreen® Dye has been published in hundreds of journal articles. This is a selection of publications using EvaGreen® Dye in a variety of applications.

EvaGreen® Dye by Application

- [Real-time qPCR](#) p. 1
- [High Resolution Melt Analysis \(HRM\)](#) p. 5
- [Digital PCR & Droplet Digital PCR \(ddPCR\)](#) p. 7
- [Fluidigm Instruments](#) p. 8
- [Isothermal Amplification](#) p. 10
- [Capillary Electrophoresis](#) p. 11
- [DNA Quantitation](#) p. 12
- [Other Applications](#) p. 12

EvaGreen® Dye References by Application

Real-time qPCR

Almeida, L. O., et al. [Accumulated SET protein up-regulates and interacts with hnRNPK, increasing its binding to nucleic acids, the Bcl-xS repression, and cellular proliferation](#). Biochem Biophys Res Commun 445(1), 196-202, (2014), DOI: 10.1016/j.bbrc.2014.01.175 (Fast EvaGreen Master Mix)

Babelova, A., et al. [Activation of Rac-1 and RhoA contributes to podocyte injury in chronic kidney disease](#). PLoS One 8(11), e80328, (2013), DOI: 10.1371/journal.pone.0080328 (Fast Plus EvaGreen Master Mix)

Camargo Lde, L., et al. [Endo-PDI is required for TNFalpha-induced angiogenesis](#). Free Radic Biol Med 65, 1398-1407, (2013), DOI: 10.1016/j.freeradbiomed.2013.09.028 (Fast Plus EvaGreen Master Mix)

Canier, L., et al. [An innovative tool for moving malaria PCR detection of parasite reservoir into the field](#). Malar J 12, 405, (2013), DOI: 10.1186/1475-2875-12-405

Chan, L. T., et al. [MicroRNAs: new insights into the pathogenesis of endodontic periapical disease](#). J Endod 39(12), 1498-1503, (2013), DOI: 10.1016/j.joen.2013.08.032 (Fast EvaGreen Master Mix)

Chen, D., et al. [High-sensitivity PCR method for detecting BRAF V600E mutations in metastatic colorectal cancer using LNA/DNA chimeras to block wild-type alleles](#). Anal Bioanal Chem, (2014), DOI: 10.1007/s00216-014-7618-x

Chen, Y., et al. [Identification of cervidae DNA in feedstuff using a real-time polymerase chain reaction method with the new fluorescence intercalating dye EvaGreen](#). J AOAC Int 92(1), 175-180, (2009),

Cheng, J., et al. [Development of a single-tube multiplex real-time PCR for detection and identification of five pathogenic targets by using melting-curve analysis with EvaGreen](#). Arch Virol, (2012), DOI: 10.1007/s00705-012-1493-6

- Choi, W. G., et al. [Salt stress-induced Ca²⁺ waves are associated with rapid, long-distance root-to-shoot signaling in plants](#). Proc Natl Acad Sci U S A 111(17), 6497-6502, (2014), DOI: 10.1073/pnas.1319955111 (Fast EvaGreen Master Mix)
- Eischeid, A. C. [SYTO dyes and EvaGreen outperform SYBR Green in real-time PCR](#). BMC Res Notes 4, 263, (2011), DOI: 10.1186/1756-0500-4-263
- Flores Bueso, Y., et al. [Characterization of FFPE-induced bacterial DNA damage and development of a repair method](#). Biol Methods Protoc 5(1), bpa015, (2020), DOI: 10.1093/biomet/bpa015
- Ganguli, A., et al. [Hands-free smartphone-based diagnostics for simultaneous detection of Zika, Chikungunya, and Dengue at point-of-care](#). Biomed Microdevices 19(4), 73, (2017), DOI: 10.1007/s10544-017-0209-9
- Gonzalez-Gonzalez, E., et al. [Portable and accurate diagnostics for COVID-19: Combined use of the miniPCR thermocycler and a well-plate reader for SARS-CoV-2 virus detection](#). PLoS One 15(8), e0237418, (2020), DOI: 10.1371/journal.pone.0237418
- Haines, A. M., et al. [Optimization of Diamond Nucleic Acid Dye for quantitative PCR](#). Biotechniques 61(4), 183-189, (2016), DOI: 000114458 [pii] 10.2144/000114458
- He, P., et al. [Multiplex real-time PCR assay for detection of pathogenic Vibrio parahaemolyticus strains](#). Mol Cell Probes, (2014), DOI: 10.1016/j.mcp.2014.06.001
- Hernandez-Arteaga, S. and Lopez-Revilla, R. [Quantitation of human papillomavirus type 16 E6 oncogene sequences by real-time or quantitative PCR with EvaGreen](#). Anal Biochem 380(1), 131-133, (2008), DOI: 10.1016/j.ab.2008.05.012
- Hu, Z., et al. [Development of a single-tube duplex EvaGreen real-time PCR for the detection and identification of EHV-1 and EHV-4](#). Appl Microbiol Biotechnol 98(9), 4179-4186, (2014), DOI: 10.1007/s00253-014-5626-6
- Khan, S. A., et al. [Detection of aacA-aphD, qacEdelta1, marA, floR, and tetA genes from multidrug-resistant bacteria: Comparative analysis of real-time multiplex PCR assays using EvaGreen\(\(R\)\) and SYBR\(\(R\)\) Green I dyes](#). Mol Cell Probes 25(2-3), 78-86, (2011), DOI: 10.1016/j.mcp.2011.01.004
- Kolm, C., et al. [DNA aptamers against bacterial cells can be efficiently selected by a SELEX process using state-of-the art qPCR and ultra-deep sequencing](#). Sci Rep 10(1), 20917, (2020), DOI: 10.1038/s41598-020-77221-9
- Kropotova, E. S., et al. [Altered expression of multiple genes involved in retinoic acid biosynthesis in human colorectal cancer](#). Pathol Oncol Res 20(3), 707-717, (2014), DOI: 10.1007/s12253-014-9751-4 (Fast Plus EvaGreen Master Mix)
- Lee, H. G., et al. [PCR-coupled Paper-based Surface-enhanced Raman Scattering \(SERS\) Sensor for Rapid and Sensitive Detection of Respiratory Bacterial DNA](#). Sensors and Actuators B: Chemical 326, (2021), DOI: 10.1016/j.snb.2020.128802
- Li, J., et al. [The Cloning and Characterization of the Enolase2 Gene of Gekko japonicus and Its Polyclonal Antibody Preparation](#). Int J Mol Sci 14(5), 8787-8800, (2013), DOI: 10.3390/ijms14058787 (Fast EvaGreen Master Mix)

Luo, Y., et al. [Massively parallel single-molecule telomere length measurement with digital real-time PCR](#). *Sci Adv* 6(34), (2020), DOI: 10.1126/sciadv.abb7944

Mao, F., et al. [Characterization of EvaGreen and the implication of its physicochemical properties for qPCR applications](#). *BMC Biotechnol* 7, 76, (2007), DOI: 10.1186/1472-6750-7-76

Moniri, A., et al. [Amplification Curve Analysis: Data-Driven Multiplexing Using Real-Time Digital PCR](#). *Anal Chem* 92(19), 13134-13143, (2020), DOI: 10.1021/acs.analchem.0c02253

Montgomery, J. L. and Wittwer, C. T. [Influence of PCR reagents on DNA polymerase extension rates measured on real-time PCR instruments](#). *Clin Chem* 60(2), 334-340, (2014), DOI: 10.1373/clinchem.2013.212829

Mossbrugger, I., et al. [EvaGreen based real-time RT-PCR assay for broad-range detection of hantaviruses in the field](#). *J Clin Virol* 58(1), 334-335, (2013), DOI: 10.1016/j.jcv.2013.06.023 qPCR to detect hantavirus

Nanayakkara, I. A., et al. [Simplifying Nucleic Acid Amplification from Whole Blood with Direct Polymerase Chain Reaction on Chitosan Microparticles](#). *Anal Chem* 89(6), 3773-3779, (2017), DOI: 10.1021/acs.analchem.7b00274

Naqvi, A. R., et al. [MicroRNAs responsive to *Aggregatibacter actinomycetemcomitans* and *Porphyromonas gingivalis* LPS modulate expression of genes regulating innate immunity in human macrophages](#). *Innate Immun* 20(5), 540-551, (2013), DOI: 10.1177/1753425913501914 (Fast EvaGreen Master Mix)

Novak, L., et al. [An integrated fluorescence detection system for lab-on-a-chip applications](#). *Lab Chip* 7(1), 27-29, (2007), DOI: 10.1039/b6111745g

Oliveira-Pelegrin, G. R., et al. [Cleaved caspase-3 expression in hypothalamic magnocellular neurons may affect vasopressin secretion during experimental polymicrobial sepsis](#). *J Neuroimmunol* 258(1-2), 10-16, (2013), DOI: 10.1016/j.jneuroim.2013.02.007 (Fast EvaGreen Master Mix)

Padirac, A., et al. [Spatial waves in synthetic biochemical networks](#). *J Am Chem Soc* 135(39), 14586-14592, (2013), DOI: 10.1021/ja403584p

Purhonen, J., et al. [A sensitive assay for dNTPs based on long synthetic oligonucleotides, EvaGreen dye and inhibitor-resistant high-fidelity DNA polymerase](#). *Nucleic Acids Res* 48(15), e87, (2020), DOI: 10.1093/nar/gkaa516

Qian, Z., et al. [Quantification of GRAF gene expression in patients with acute myeloid leukemia using EvaGreen real time quantitative PCR](#). *Zhonghua Yi Xue Yi Chuan Xue Za Zhi* 27(3), 290-293, (2010), DOI: 10.3760/cma.j.issn.1003-9406.2010.0.012

Rodrigues, E. F., et al. [Epigenetic alterations of p15\(INK4B\) and p16\(INK4A\) genes in pediatric primary myelodysplastic syndrome](#). *Leuk Lymphoma* 51(10), 1887-1894, (2010), DOI: 10.3109/10428194.2010.505820 (Fast EvaGreen Master Mix)

Sang, F., et al. [Development of a high-throughput real time PCR based on a hot-start alternative for Pfu mediated by quantum dots](#). *Nanoscale* 7(38), 15852-15862, (2015), DOI: 10.1039/c5nr03596a

- Sang, F., et al. [CdTe quantum dots enhance feasibility of EvaGreen-based real-time PCR with decent amplification fidelity](#). Mol Biotechnol 54(3), 969-976, (2013), DOI: 10.1007/s12033-013-9650-z
- Shcheglovitov, A., et al. [SHANK3 and IGF1 restore synaptic deficits in neurons from 22q13 deletion syndrome patients](#). Nature 503(7475), 267-271, (2013), DOI: 10.1038/nature12618
- Shi, W., et al. [EvaGreen-based real-time PCR assay for sensitive detection of salmonid alphavirus](#). Mol Cell Probes 39, 7-13, (2018), DOI: 10.1016/j.mcp.2018.03.002
- Shu, B., et al. [A sample-to-answer, real-time convective polymerase chain reaction system for point-of-care diagnostics](#). Biosens Bioelectron 97, 360-368, (2017), DOI: 10.1016/j.bios.2017.06.014
- Swennenhuis, J. F., et al. [Efficiency of whole genome amplification of single circulating tumor cells enriched by CellSearch and sorted by FACS](#). Genome Med 5(11), 106, (2013), DOI: 10.1186/gm510
- Timon-Gomez, A., et al. [Differential regulation of mitochondrial pyruvate carrier genes modulates respiratory capacity and stress tolerance in yeast](#). PLoS One 8(11), e79405, (2013), DOI: 10.1371/journal.pone.0079405 (Fast EvaGreen Master Mix)
- Velez, D. O., et al. [Massively parallel digital high resolution melt for rapid and absolutely quantitative sequence profiling](#). Sci Rep 7, 42326, (2017), DOI: 10.1038/srep42326
- Veltkamp, H. W., et al. [Disposable DNA Amplification Chips with Integrated Low-Cost Heaters](#). Micromachines (Basel) 11(3), (2020), DOI: 10.3390/mi11030238
- Vig, P. J., et al. [Knockdown of acid-sensing ion channel 1a \(ASIC1a\) suppresses disease phenotype in SCA1 mouse model](#). Cerebellum 13(4), 479-490, (2014), DOI: 10.1007/s12311-014-0563-6 (Fast EvaGreen Master Mix)
- Wang, C., et al. [Rapid screening of MMACHC gene mutations by high-resolution melting curve analysis](#). Mol Genet Genomic Med 8(6), e1221, (2020), DOI: 10.1002/mgg3.1221
- Wang, C. P., et al. [Isoquercetin protects cortical neurons from oxygen-glucose deprivation-reperfusion induced injury via suppression of TLR4-NF-small ka, CyrillicB signal pathway](#). Neurochem Int 63(8), 741-749, (2013), DOI: 10.1016/j.neuint.2013.09.018 (Fast Plus EvaGreen Master Mix)
- Wei, C., et al. [Identification of the crucial molecular events during the large-scale myoblast fusion in sheep](#). Physiol Genomics 46(12), 429-440, (2014), DOI: 10.1152/physiolgenomics.00184.2013 (Fast EvaGreen Master Mix)
- Zhang, H., et al. [PCR Multiplexing Based on a Single Fluorescent Channel Using Dynamic Melting Curve Analysis](#). ACS Omega 5(46), 30267-30273, (2020), DOI: 10.1021/acsomega.0c04766
- Zhang, Q., et al. [Involvement of ERK1/2 pathway in neuroprotective effects of pyrroloquinoline quinine against rotenone-induced SH-SY5Y cell injury](#). Neuroscience 270, 183-191, (2014), DOI: 10.1016/j.neuroscience.2014.04.022 (Fast EvaGreen Master Mix)
- Zhu, S. and Coffman, J. A. [Simple and fast quantification of DNA damage by real-time PCR, and its application to nuclear and mitochondrial DNA from multiple tissues of aging zebrafish](#). BMC Res Notes 10(1), 269, (2017), DOI: 10.1186/s13104-017-2593-x

High Resolution Melt Analysis (HRM)

Athamanolap, P., et al. [Droplet Array Platform for High-Resolution Melt Analysis of DNA Methylation Density](#). J Lab Autom, (2013), DOI: 10.1177/2211068213507923

Cai, X. Q., et al. [Rapid detection and simultaneous genotyping of Cronobacter spp. \(formerly Enterobacter sakazakii\) in powdered infant formula using real-time PCR and high resolution melting \(HRM\) analysis](#). PLoS One 8(6), e67082, (2013), DOI: 10.1371/journal.pone.0067082 (Fast Plus EvaGreen Master Mix)

Cubas-Atienzar, A. I., et al. [A novel air-dried multiplex high resolution melt assay for detection of extended spectrum beta-lactamase and carbapenemase genes](#). MedRxiv, (2021), DOI: 10.1101/2021.02.18.21251950

Dhami, M. K. and Kumarasinghe, L. [A HRM real-time PCR assay for rapid and specific identification of the emerging pest spotted-wing drosophila \(Drosophila suzukii\)](#). PLoS One 9(6), e98934, (2014), DOI: 10.1371/journal.pone.0098934

Dong, J., et al. [Single-color multiplexing by the integration of high-resolution melting pattern recognition with loop-mediated isothermal amplification](#). Chem Commun (Camb) 55(17), 2457-2460, (2019), DOI: 10.1039/c8cc09741k

Flores Bueso, Y., et al. [Characterization of FFPE-induced bacterial DNA damage and development of a repair method](#). Biol Methods Protoc 5(1), bpaa015, (2020), DOI: 10.1093/biomethods/bpaa015

Fraley, S. I., et al. [Universal digital high-resolution melt: a novel approach to broad-based profiling of heterogeneous biological samples](#). Nucleic Acids Res 41(18), e175, (2013), DOI: 10.1093/nar/gkt684 U

Hanson, E. K. and Ballantyne, J. [Rapid and inexpensive body fluid identification by RNA profiling-based multiplex High Resolution Melt \(HRM\) analysis](#). F1000Res 2, 281, (2013), DOI: 10.12688/f1000research.2-281.v2

Kozina, V., et al. [A one-step real-time multiplex PCR for screening Y-chromosomal microdeletions without downstream amplicon size analysis](#). PLoS One 6(8), e23174, (2011), DOI: 10.1371/journal.pone.0023174

Li, Y. D., et al. [A cost-effective high-resolution melting approach using the EvaGreen dye for DNA polymorphism detection and genotyping in plants](#). J Integr Plant Biol 52(12), 1036-1042, (2010), DOI: 10.1111/j.1744-7909.2010.01001.x

Liyanage, K. E., et al. [High-resolution melting analysis for detection of familial ligand-defective apolipoprotein B-100 mutations](#). Ann Clin Biochem 45(Pt 2), 170-176, (2008), DOI: 10.1258/acb.2007.007077

Mader, E., et al. [Quantitative high-resolution melting analysis for detecting adulterations](#). Anal Biochem 409(1), 153-155, (2011), DOI: 10.1016/j.ab.2010.10.009

Maltese, P., et al. [High resolution melting \(HRM\) analysis for the detection of ER22/23EK, BclI, and N363S polymorphisms of the glucocorticoid receptor gene](#). J Steroid Biochem Mol Biol 113(3-5), 269-274, (2009), DOI: 10.1016/j.jsbmb.2009.01.012

- Miller, M., et al. [High-Resolution Melting Curve Analysis for Identification of Pasteurellaceae Species in Experimental Animal Facilities](#). PLoS One 10(11), e0142560, (2015), DOI: 10.1371/journal.pone.0142560 PONE-D-15-35081 [pii]
- Nemcova, E., et al. [Rapid identification of medically important Candida isolates using high resolution melting analysis](#). PLoS One 10(2), e0116940, (2015), DOI: 10.1371/journal.pone.0116940 PONE-D-14-41648 [pii]
- Pisanic, T. R., 2nd, et al. [DREAMing: a simple and ultrasensitive method for assessing intratumor epigenetic heterogeneity directly from liquid biopsies](#). Nucleic Acids Res 43(22), e154, (2015), DOI: gkv795 [pii] 10.1093/nar/gkv795
- Radvanszky, J., et al. [Comparison of different DNA binding fluorescent dyes for applications of high-resolution melting analysis](#). Clin Biochem 48(9), 609-616, (2015), DOI: S0009-9120(15)00032-6 [pii] 10.1016/j.clinbiochem.2015.01.010
- Rao, P., et al. [Development of an EvaGreen-based multiplex real-time PCR assay with melting curve analysis for simultaneous detection and differentiation of six viral pathogens of porcine reproductive and respiratory disorder](#). J Virol Methods 208, 56-62, (2014), DOI: 10.1016/j.jviromet.2014.06.027
- Ronai, Z., et al. [Molecular identification of Mycobacterium avium subsp. silvaticum by duplex high-resolution melt analysis and subspecies-specific real-time PCR](#). J Clin Microbiol 53(5), 1582-1587, (2015), DOI: JCM.03556-14 [pii] 10.1128/JCM.03556-14
- Tanaka, M., et al. [High-resolution melting analysis for genotyping Duffy, Kidd and Diego blood group antigens](#). Leg Med (Tokyo) 13(1), 1-6, (2011), DOI: 10.1016/j.legalmed.2010.08.004
- Twist, G. P., et al. [High-resolution melt analysis to detect sequence variations in highly homologous gene regions: application to CYP2B6](#). Pharmacogenomics 14(8), 913-922, (2013), DOI: 10.2217/pgs.13.66
- Valenzuela-Munoz, V., et al. [SNP discovery and High Resolution Melting Analysis from massive transcriptome sequencing in the California red abalone Haliotis rufescens](#). Mar Genomics 10, 11-16, (2013), DOI: 10.1016/j.margen.2012.12.003 (Fast EvaGreen Master Mix)
- Velez, D. O., et al. [Massively parallel digital high resolution melt for rapid and absolutely quantitative sequence profiling](#). Sci Rep 7, 42326, (2017), DOI: 10.1038/srep42326
- Wang, C., et al. [Rapid screening of MMACHC gene mutations by high-resolution melting curve analysis](#). Mol Genet Genomic Med 8(6), e1221, (2020), DOI: 10.1002/mgg3.1221
- Wang, J., et al. [Assessment for Melting Temperature Measurement of Nucleic Acid by HRM](#). J Anal Methods Chem 2016, 5318935, (2016), DOI: 10.1155/2016/5318935
- White, H. E., et al. [Methylation-sensitive high-resolution melting-curve analysis of the SNRPN gene as a diagnostic screen for Prader-Willi and Angelman syndromes](#). Clin Chem 53(11), 1960-1962, (2007), DOI: 10.1373/clinchem.2007.093351
- Xu, X., et al. [Rapid screening of UPB1 gene variations by high resolution melting curve analysis](#). Exp Ther Med 21(4), 403, (2021), DOI: 10.3892/etm.2021.9834

You, C. G., et al. [Association analysis of single nucleotide polymorphisms of proinflammatory cytokine and their receptors genes with rheumatoid arthritis in northwest Chinese Han population](#). Cytokine 61(1), 133-138, (2013), DOI: 10.1016/j.cyto.2012.09.007 (Fast EvaGreen Master Mix)

Digital PCR & Droplet Digital PCR (ddPCR)

Athamanolap, P., et al. [Droplet Array Platform for High-Resolution Melt Analysis of DNA Methylation Density](#). J Lab Autom, (2013), DOI: 10.1177/2211068213507923

Atour, A., et al. [iSpinach: a fluorogenic RNA aptamer optimized for in vitro applications](#). Nucleic Acids Res 44(6), 2491-2500, (2016), DOI: gkw083 [pii] 10.1093/nar/gkw083

Falzone, L., et al. [Sensitivity assessment of droplet digital PCR for SARS-CoV-2 detection](#). Int J Mol Med 46(3), 957-964, (2020), DOI: 10.3892/ijmm.2020.4673

Fraley, S. I., et al. [Universal digital high-resolution melt: a novel approach to broad-based profiling of heterogeneous biological samples](#). Nucleic Acids Res 41(18), e175, (2013), DOI: 10.1093/nar/gkt684

McDermott, G. P., et al. [Multiplexed target detection using DNA-binding dye chemistry in droplet digital PCR](#). Anal Chem 85(23), 11619-11627, (2013), DOI: 10.1021/ac403061n

Miotke, L. K., et al. [High Sensitivity Detection and Quantitation of DNA Copy Number and Single Nucleotide Variants with Single Color Droplet Digital PCR](#). Anal Chem, (2014), DOI: 10.1021/ac403843j

Moniri, A., et al. [High-Level Multiplexing in Digital PCR with Intercalating Dyes by Coupling Real-Time Kinetics and Melting Curve Analysis](#). Anal Chem 92(20), 14181-14188, (2020), DOI: 10.1021/acs.analchem.0c03298

Moniri, A., et al. [Amplification Curve Analysis: Data-Driven Multiplexing Using Real-Time Digital PCR](#). Anal Chem 92(19), 13134-13143, (2020), DOI: 10.1021/acs.analchem.0c02253

Ryckelynck, M., et al. [Using droplet-based microfluidics to improve the catalytic properties of RNA under multiple-turnover conditions](#). RNA 21(3), 458-469, (2015), DOI:10.1261/rna.048033.114

Sidore, A. M., et al. [Enhanced sequencing coverage with digital droplet multiple displacement amplification](#). Nucleic Acids Res 44(7), e66, (2016), DOI: 10.1093/nar/gkv1493

Wang, Y., et al. [Digital PCR using micropatterned superporous absorbent array chips](#). Analyst 141(12), 3821-3831, (2016), DOI: 10.1039/c6an00164e

Wood-Bouwens, C., et al. [Single-Color Digital PCR Provides High-Performance Detection of Cancer Mutations from Circulating DNA](#). J Mol Diagn 19(5), 697-710, (2017), DOI: 10.1016/j.jmoldx.2017.05.003

Zhang, H., et al. [Isolation and Analysis of Rare Norovirus Recombinants from Coinfected Mice Using Drop-Based Microfluidics](#). J Virol 89(15), 7722-7734, (2015), DOI: 10.1128/JVI.01137-15

Zhang, H., et al. [PCR Multiplexing Based on a Single Fluorescent Channel Using Dynamic Melting Curve Analysis](#). ACS Omega 5(46), 30267-30273, (2020), DOI: 10.1021/acsomega.0c04766

Fluidigm Instruments

Adamski, M. G., et al. [Next-generation qPCR for the high-throughput measurement of gene expression in multiple leukocyte subsets](#). J Biomol Screen 18(9), 1008-1017, (2013), DOI: 10.1177/1087057113489882 (Fast EvaGreen Master Mix)

Allen, E. K., et al. [Evaluation of replication of variants associated with genetic risk of otitis media](#). PLoS One 9(8), e104212, (2014), DOI: 10.1371/journal.pone.0104212 (Fast Probe Master Mix)

Barribeau, S. M. and Schmid-Hempel, P. [Qualitatively different immune response of the bumblebee host, *Bombus terrestris*, to infection by different genotypes of the trypanosome gut parasite, *Crithidia bombi*](#). Infect Genet Evol 20, 249-256, (2013), DOI: 10.1016/j.meegid.2013.09.014

Brunner, F. S., et al. [Immune gene expression in *Bombus terrestris*: signatures of infection despite strong variation among populations, colonies, and sister workers](#). PLoS One 8(7), e68181, (2013), DOI: 10.1371/journal.pone.0068181

Bugarel, M., et al. [Molecular identification in monophasic and nonmotile variants of *Salmonella enterica* serovar Typhimurium](#). Microbiologyopen 1(4), 481-489, (2012), DOI: 10.1002/mbo3.39

Citri, A., et al. [Comprehensive qPCR profiling of gene expression in single neuronal cells](#). Nat Protoc 7(1), 118-127, (2012), DOI: 10.1038/nprot.2011.430

Diepeveen, E. T., et al. [Immune-related functions of the *Hivep* gene family in East African cichlid fishes](#). G3 (Bethesda) 3(12), 2205-2217, (2013), DOI: 10.1534/g3.113.008839

Flynn, J. M., et al. [Single cell gene expression profiling of cortical osteoblast lineage cells](#). Bone 53(1), 174-181, (2013), DOI: 10.1016/j.bone.2012.11.043

Guo, G., et al. [Mapping cellular hierarchy by single-cell analysis of the cell surface repertoire](#). Cell Stem Cell 13(4), 492-505, (2013), DOI: 10.1016/j.stem.2013.07.017

Ho, D. W., et al. [Gene expression profiling of liver cancer stem cells by RNA-sequencing](#). PLoS One 7(5), e37159, (2012), DOI: 10.1371/journal.pone.0037159

Isom, S. C., et al. [Transcriptional profiling by RNA-Seq of peri-attachment porcine embryos generated by a variety of assisted reproductive technologies](#). Physiol Genomics 45(14), 577-589, (2013), DOI: 10.1152/physiolgenomics.00094.2012

Jasnos, L., et al. [Identifying division symmetry of mouse embryonic stem cells: negative impact of DNA methyltransferases on symmetric self-renewal](#). Stem Cell Reports 1(4), 360-369, (2013), DOI: 10.1016/j.stemcr.2013.08.005

Jasnos, L. and Sawado, T. [Determining cell division symmetry through the dissection of dividing cells using single-cell expression analysis](#). Nat Protoc 9(3), 505-516, (2014), DOI: 10.1038/nprot.2014.032

Johnston, W. T., et al. [Relationship between *Plasmodium falciparum* malaria prevalence, genetic diversity and endemic Burkitt lymphoma in Malawi](#). Sci Rep 4, 3741, (2014), DOI: 10.1038/srep03741 (Fast Probe Master Mix)

Kerenyi, M. A., et al. [Histone demethylase *Lsd1* represses hematopoietic stem and progenitor cell signatures during blood cell maturation](#). Elife 2, e00633, (2013), DOI: 10.7554/eLife.00633

- Khan, S. A., et al. [The mQTL hotspot on linkage group 16 for phenolic compounds in apple fruits is probably the result of a leucoanthocyanidin reductase gene at that locus](#). BMC Res Notes 5, 618, (2012), DOI: 10.1186/1756-0500-5-618
- Laurell, H., et al. [Correction of RT-qPCR data for genomic DNA-derived signals with ValidPrime](#). Nucleic Acids Res 40(7), e51, (2012), DOI: 10.1093/nar/gkr1259
- Liang-Chu, M. M., et al. [Human biosample authentication using the high-throughput, cost-effective SNPtrace\(TM\) system](#). PLoS One 10(2), e0116218, (2015), DOI: 10.1371/journal.pone.0116218 PONE-D-14-45036 [pii]
- May-Panloup, P., et al. [Molecular characterization of corona radiata cells from patients with diminished ovarian reserve using microarray and microfluidic-based gene expression profiling](#). Hum Reprod 27(3), 829-843, (2012), DOI: 10.1093/humrep/der431
- Murima, P., et al. [Exploring the mode of action of bioactive compounds by microfluidic transcriptional profiling in mycobacteria](#). PLoS One 8(7), e69191, (2013), DOI: 10.1371/journal.pone.0069191
- Pasca, S. P., et al. [Using iPSC-derived neurons to uncover cellular phenotypes associated with Timothy syndrome](#). Nat Med 17(12), 1657-1662, (2011), DOI: 10.1038/nm.2576
- Pedersen, R., et al. [Characterisation of gut microbiota in Ossabaw and Gottingen minipigs as models of obesity and metabolic syndrome](#). PLoS One 8(2), e56612, (2013), DOI: 10.1371/journal.pone.0056612
- Piazza, R. M., et al. [Molecular and phenotypic characterization of Escherichia coli O26:H8 among diarrheagenic E. coli O26 strains isolated in Brazil](#). Appl Environ Microbiol 79(22), 6847-6854, (2013), DOI: 10.1128/AEM.01693-13
- Rodgaard, T., et al. [Cloning changes the response to obesity of innate immune factors in blood, liver, and adipose tissues in domestic pigs](#). Cell Reprogram 15(3), 185-194, (2013), DOI: 10.1089/cell.2012.0091
- Rusnakova, V., et al. [Heterogeneity of astrocytes: from development to injury - single cell gene expression](#). PLoS One 8(8), e69734, (2013), DOI: 10.1371/journal.pone.0069734
- Schunter, C., et al. [SNP development from RNA-seq data in a nonmodel fish: how many individuals are needed for accurate allele frequency prediction?](#). Mol Ecol Resour 14(1), 157-165, (2014), DOI: 10.1111/1755-0998.12155 (Fast Probe Master Mix)
- Shcheglovitov, A., et al. [SHANK3 and IGF1 restore synaptic deficits in neurons from 22q13 deletion syndrome patients](#). Nature 503(7475), 267-271, (2013), DOI: 10.1038/nature12618
- Skovgaard, K., et al. [Expression of innate immune genes, proteins and microRNAs in lung tissue of pigs infected experimentally with influenza virus \(H1N2\)](#). Innate Immun 19(5), 531-544, (2013), DOI: 10.1177/1753425912473668
- Sreenivasan, R., et al. [Gonad Differentiation in Zebrafish Is Regulated by the Canonical Wnt Signaling Pathway](#). Biol Reprod, (2013), DOI: 10.1095/biolreprod.113.110874
- Veltkamp, H. W., et al. [Disposable DNA Amplification Chips with Integrated Low-Cost Heaters](#). Micromachines (Basel) 11(3), (2020), DOI: 10.3390/mi11030238

Weinstein, J. A., et al. [Correlation of gene expression and genome mutation in single B-cells](#). PLoS One 8(6), e67624, (2013), DOI: 10.1371/journal.pone.0067624

Wopereis, S., et al. [Assessment of inflammatory resilience in healthy subjects using dietary lipid and glucose challenges](#). BMC Med Genomics 6, 44, (2013), DOI: 10.1186/1755-8794-6-44

Isothermal Amplification

Alekseenko, A., et al. [Direct detection of SARS-CoV-2 using non-commercial RT-LAMP reagents on heat-inactivated samples](#). Sci Rep 11(1), 1820, (2021), DOI: 10.1038/s41598-020-80352-8

Chen, W., et al. [Mobile Platform for Multiplexed Detection and Differentiation of Disease-Specific Nucleic Acid Sequences, Using Microfluidic Loop-Mediated Isothermal Amplification and Smartphone Detection](#). Anal Chem 89(21), 11219-11226, (2017), DOI: 10.1021/acs.analchem.7b02478

Dong, J., et al. [Single-color multiplexing by the integration of high-resolution melting pattern recognition with loop-mediated isothermal amplification](#). Chem Commun (Camb) 55(17), 2457-2460, (2019), DOI: 10.1039/c8cc09741k

Duarte, C., et al. [On-chip parallel detection of foodborne pathogens using loop-mediated isothermal amplification](#). Biomed Microdevices, (2013), DOI: 10.1007/s10544-013-9769-5

El-Tholoth, M., et al. [A Single and Two-Stage, Closed-Tube, Molecular Test for the 2019 Novel Coronavirus \(COVID-19\) at Home, Clinic, and Points of Entry](#). ChemRxiv, (2020), DOI: 10.26434/chemrxiv.11860137.v1 10.26434/chemrxiv.11860137

Fukuta, S., et al. [Development of loop-mediated isothermal amplification assay for the detection of Pythium myriotylum](#). Lett Appl Microbiol 59(1), 49-57, (2014), DOI: 10.1111/lam.12244

Ganguli, A., et al. [Rapid isothermal amplification and portable detection system for SARS-CoV-2](#). Proc Natl Acad Sci U S A 117(37), 22727-22735, (2020), DOI: 10.1073/pnas.2014739117

Ganguli, A., et al. [Reverse Transcription Loop-Mediated Isothermal Amplification Assay for Ultrasensitive Detection of SARS-CoV-2 in Saliva and Viral Transport Medium Clinical Samples](#). Anal Chem 93(22), 7797-7807, (2021), DOI: 10.1021/acs.analchem.0c05170

Ganguli, A., et al. [Hands-free smartphone-based diagnostics for simultaneous detection of Zika, Chikungunya, and Dengue at point-of-care](#). Biomed Microdevices 19(4), 73, (2017), DOI: 10.1007/s10544-017-0209-9

Garcia, D. C., et al. [Expression of urokinase type plasminogen activator receptor \(uPAR\) in the bovine oviduct: Relationship with uPA effect on oviductal epithelial cells](#). Res Vet Sci 97(1), 118-123, (2014), DOI: 10.1016/j.rvsc.2014.05.012 (Fast Plus EvaGreen Master Mix)

Goldmeyer, J., et al. [Development of a novel one-tube isothermal reverse transcription thermophilic helicase-dependent amplification platform for rapid RNA detection](#). J Mol Diagn 9(5), 639-644, (2007), DOI: 10.2353/jmoldx.2007.070012

Li, Y., et al. [Primase-based whole genome amplification](#). Nucleic Acids Res 36(13), e79, (2008), DOI: 10.1093/nar/gkn377

- Liao, S. C., et al. [Smart Cup: A Minimally-Instrumented, Smartphone-Based Point-of-Care Molecular Diagnostic Device](#). *Sens Actuators B Chem* 229, 232-238, (2016), DOI: 10.1016/j.snb.2016.01.073
- Liu, C., et al. [Membrane-based, sedimentation-assisted plasma separator for point-of-care applications](#). *Anal Chem* 85(21), 10463-10470, (2013), DOI: 10.1021/ac402459h
- Loo, J. F. C., et al. [An aptamer-based bio-barcode assay with isothermal recombinase polymerase amplification for cytochrome-c detection and anti-cancer drug screening](#). *Talanta* 115(0), 159-165, (2013), DOI: 10.1016/j.talanta.2013.04.051
- Montgomery, J. L., et al. [Stopped-flow DNA polymerase assay by continuous monitoring of dNTP incorporation by fluorescence](#). *Anal Biochem* 441(2), 133-139, (2013), DOI: 10.1016/j.ab.2013.07.008
- Salm, E., et al. [Electrical detection of nucleic acid amplification using an on-chip quasi-reference electrode and a PVC REFET](#). *Anal Chem* 86(14), 6968-6975, (2014), DOI: 10.1021/ac500897t
- Schaerli, Y., et al. [Isothermal DNA amplification using the T4 replisome: circular nicking endonuclease-dependent amplification and primase-based whole-genome amplification](#). *Nucleic Acids Res* 38(22), e201, (2010), DOI: 10.1093/nar/gkq795
- Wang, D. and Liu, Y. [Development of Primer Sets for Loop-Mediated Isothermal Amplification that Enables Rapid and Specific Detection of *Streptococcus dysgalactiae*, *Streptococcus uberis* and *Streptococcus agalactiae*](#). *Int J Environ Res Public Health* 12(6), 5735-5742, (2015), DOI: ijerph120605735 [pii] 10.3390/ijerph120605735
- Wang, D., et al. [A Comparison of In-House Real-Time LAMP Assays with a Commercial Assay for the Detection of Pathogenic Bacteria](#). *Molecules* 20(6), 9487-9495, (2015), DOI: molecules20069487 [pii] 10.3390/molecules20069487
- Wang, D. G., et al. [Two methods for increased specificity and sensitivity in loop-mediated isothermal amplification](#). *Molecules* 20(4), 6048-6059, (2015), DOI: molecules20046048 [pii] 10.3390/molecules20046048
- Yi, C., et al. [Nanoporous hydrogel for direct digital nucleic acid amplification in untreated complex matrices for single bacteria counting](#). *Biosens Bioelectron* 184, 113199, (2021), DOI: 10.1016/j.bios.2021.113199
- Zhou, Q. J., et al. [Development and evaluation of a real-time fluorogenic loop-mediated isothermal amplification assay integrated on a microfluidic disc chip \(on-chip LAMP\) for rapid and simultaneous detection of ten pathogenic bacteria in aquatic animals](#). *J Microbiol Methods* 104, 26-35, (2014), DOI: 10.1016/j.mimet.2014.06.008

Capillary Electrophoresis

- Sang, F., et al. [Genetic mutation analysis by CE with LIF detection using inverse-flow derivatization of DNA fragments](#). *Electrophoresis* 27(19), 3846-3855, (2006), DOI: 10.1002/elps.200600160
- Sang, F. and Ren, J. [Capillary electrophoresis of double-stranded DNA fragments using a new fluorescence intercalating dye EvaGreen](#). *J Sep Sci* 29(9), 1275-1280, (2006), DOI: 10.1002/jssc.200600029

DNA Quantification

Ihrig, J., et al. [Application of the DNA-specific dye EvaGreen for the routine quantification of DNA in microplates](#). *Anal Biochem* 359(2), 265-267, (2006), DOI: 10.1016/j.ab.2006.07.043

Mao, F., et al. [Characterization of EvaGreen and the implication of its physicochemical properties for qPCR applications](#). *BMC Biotechnol* 7, 76, (2007), DOI: 10.1186/1472-6750-7-76

Purhonen, J., et al. [A sensitive assay for dNTPs based on long synthetic oligonucleotides, EvaGreen dye and inhibitor-resistant high-fidelity DNA polymerase](#). *Nucleic Acids Res* 48(15), e87, (2020), DOI: 10.1093/nar/gkaa516

Wang, W., et al. [DNA quantification using EvaGreen and a real-time PCR instrument](#). *Anal Biochem* 356(2), 303-305, (2006), DOI: 10.1016/j.ab.2006.05.027

Other Applications

Ahrberg, C. D., et al. [Palm-Sized Device for Point-of-Care Ebola Detection](#). *Anal Chem* 88(9), 4803-4807, (2016), DOI: 10.1021/acs.analchem.6b00278

Chiaraviglio, L. and Kirby, J. E. [Evaluation of impermeant, DNA-binding dye fluorescence as a real-time readout of eukaryotic cell toxicity in a high throughput screening format](#). *Assay Drug Dev Technol* 12(4), 219-228, (2014), DOI: 10.1089/adt.2014.577

Everitt, M. L., et al. [Sample-to-Answer Diagnostic System for the Detection of Circulating Histones in Whole Blood](#). *ACS Sens*, (2021), DOI: 10.1021/acssensors.1c00886

Gilat, N., et al. [Rapid Quantification of Oxidation and UV Induced DNA Damage by Repair Assisted Damage Detection-\(Rapid RADD\)](#). *Anal Chem* 92(14), 9887-9894, (2020), DOI: 10.1021/acs.analchem.0c01393

Huang, G., et al. [fM to aM nucleic acid amplification for molecular diagnostics in a non-stick-coated metal microfluidic bioreactor](#). *Sci Rep* 4, 7344, (2014), DOI: srep07344 [pii] 10.1038/srep07344

Luo, Y., et al. [Massively parallel single-molecule telomere length measurement with digital real-time PCR](#). *Sci Adv* 6(34), (2020), DOI: 10.1126/sciadv.abb7944

Miotke, L., et al. [Enzyme-Free Detection of Mutations in Cancer DNA Using Synthetic Oligonucleotide Probes and Fluorescence Microscopy](#). *PLoS One* 10(8), e0136720, (2015), DOI: 10.1371/journal.pone.0136720 PONE-D-15-21710 [pii]

Moyal, L., et al. [AN-7, a butyric acid prodrug, sensitizes cutaneous T-cell lymphoma cell lines to doxorubicin via inhibition of DNA double strand breaks repair](#). *Invest New Drugs* 36(1), 1-9, (2018), DOI: 10.1007/s10637-017-0500-x

Padirac, A., et al. [Spatial waves in synthetic biochemical networks](#). *J Am Chem Soc* 135(39), 14586-14592, (2013), DOI: 10.1021/ja403584p

Shoute, L. C. T. and Loppnow, G. R. [Characterization of the binding interactions between EvaGreen dye and dsDNA](#). *Phys Chem Chem Phys* 20(7), 4772-4780, (2018), DOI: 10.1039/c7cp06058k

Shu, B., et al. [A handheld flow genetic analysis system \(FGAS\): towards rapid, sensitive, quantitative and multiplex molecular diagnosis at the point-of-care level](#). Lab Chip 15(12), 2597-2605, (2015), DOI: 10.1039/c5lc00139k

Zadorin, A. S., et al. [Synthesis of programmable reaction-diffusion fronts using DNA catalyzers](#). Phys Rev Lett 114(6), 068301, (2015), DOI: 10.1103/PhysRevLett.114.068301

Zhu, G., et al. [Building Fluorescent DNA Nanodevices on Target Living Cell Surfaces](#). Angew Chem Int Ed Engl 52(21), 5490-5496, (2013), DOI: 10.1002/anie.201301439