

---

## Selected CellBrite® and MemBrite® References

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

### CellBrite® and MemBrite® References, by Product

- [CellBrite® & CellBrite® NIR Cytoplasmic Membrane Dyes](#) p. 1
- [CellBrite® Fix Membrane Stains](#) p. 4
- [MemBrite® Fix Cell Surface Staining Kits](#) p. 5

### CellBrite® and MemBrite® References for Specialized Applications

- [Organoids, Spheroids, or Matrigel/Hydrogel Cultures](#) p. 6
- [Exosome and EV Labeling](#) p. 6
- [Super Resolution Imaging](#) p. 6

## CellBrite® and MemBrite® References by Product

### CellBrite® & CellBrite® NIR Cytoplasmic Membrane Dyes

Alphonse, R. S., et al. [Existence, functional impairment, and lung repair potential of endothelial colony-forming cells in oxygen-induced arrested alveolar growth](#). *Circulation* 129(21), 2144-2157, (2014), DOI: 10.1161/CIRCULATIONAHA.114.009124 (CellBrite® Red)

Anguiano, M., et al. [The use of mixed collagen-Matrigel matrices of increasing complexity recapitulates the biphasic role of cell adhesion in cancer cell migration: ECM sensing, remodeling and forces at the leading edge of cancer invasion](#). *PLoS One* 15(1), e0220019, (2020), DOI: 10.1371/journal.pone.0220019 (CellBrite® Red, cell labeling before Matrigel® culture)

Aoyama, T., et al. [cAMP Signaling Pathway Prevents Dasatinib-Induced Vascular Hyperpermeability](#). *Biol Pharm Bull* 44(8), 1101-1110, (2021), DOI: 10.1248/bpb.b21-00270 (CellBrite® Orange)

Belmont-Monroy, L., et al. [Characterization of a novel AraC/XylS-regulated family of N-acyltransferases in pathogens of the order Enterobacterales](#). *PLoS Pathog* 16(8), e1008776, (2020), DOI: 10.1371/journal.ppat.1008776 (CellBrite® Green)

Bergenheim, F., et al. [Fluorescence-based tracing of transplanted intestinal epithelial cells using confocal laser endomicroscopy](#). *Stem Cell Res Ther* 10(1), 148, (2019), DOI: 10.1186/s13287-019-1246-5 (CellBrite® Green)

Dorrello, N. V., et al. [Functional vascularized lung grafts for lung bioengineering](#). *Sci Adv* 3(8), e1700521, (2017), DOI: 10.1126/sciadv.1700521 (CellBrite® NIR790)

Dorst, N., et al. [Analysis of cellular composition of co-culture spheroids](#). *Ann Anat* 196(5), 303-311, (2014), DOI: 10.1016/j.aanat.2014.05.038 (CellBrite® Blue)

Dryden, G. W., et al. [Comparison of Stromal Vascular Fraction with or Without a Novel Bioscaffold to Fibrin Glue in a Porcine Model of Mechanically Induced Anorectal Fistula](#). *Inflamm Bowel Dis* 23(11), 1962-1971, (2017), DOI: 10.1097/MIB.0000000000001254 (CellBrite® Red)

Garcia-Contreras, M. and Thakor, A. S. [Human adipose tissue-derived mesenchymal stem cells and their extracellular vesicles modulate lipopolysaccharide activated human microglia](#). Cell Death Discov 7(1), 98, (2021), DOI: 10.1038/s41420-021-00471-7 (CellBrite® Green, CellBrite® Orange)

Garcia-Garcia, L., et al. [Long-Term PEDF Release in Rat Iris and Retinal Epithelial Cells after Sleeping Beauty Transposon-Mediated Gene Delivery](#). Mol Ther Nucleic Acids 9, 1-11, (2017), DOI: 10.1016/j.omtn.2017.08.001 (CellBrite® Red)

Guimarães, C. F., et al. [3D flow-focusing microfluidic biofabrication: One-chip-fits-all hydrogel fiber architectures](#). Applied Materials Today 23, (2021), DOI: 10.1016/j.apmt.2021.101013 (CellBrite® Green, CellBrite® Orange, cell labeling before hydrogel culture)

Hart, M., et al. [MARK2/Par1b present at retraction fibres corrects spindle off-centering induced by actin disassembly](#). bioRxiv, (2018), DOI: 10.1101/508861 (CellBrite® Red)

Hori, T., et al. [Three-Dimensional Cell Sheet Construction Method with a Polyester Micromesh Sheet](#). Tissue Eng Part C Methods 26(3), 170-179, (2020), DOI: 10.1089/ten.tec.2019.0330 (CellBrite® Green, CellBrite® Red)

Juang, Y. P., et al. [Synthesis, distribution analysis and mechanism studies of N-acyl glucosamine-bearing oleanolic saponins](#). Bioorg Chem 99, 103835, (2020), DOI: 10.1016/j.bioorg.2020.103835 (CellBrite® Blue)

Lazaro-Ibanez, E., et al. [Metastatic state of parent cells influences the uptake and functionality of prostate cancer cell-derived extracellular vesicles](#). J Extracell Vesicles 6(1), 1354645, (2017), DOI: 10.1080/20013078.2017.1354645 (CellBrite® Green)

Li, K., et al. [Ferumoxylol-based Dual-modality Imaging Probe for Detection of Stem Cell Transplant Rejection](#). Nanotheranostics 2(4), 306-319, (2018), DOI: 10.7150/ntno.26389 (CellBrite® Red)

Luangtrakul, W., et al. [Cytotoxicity of Vibrio parahaemolyticus AHPND toxin on shrimp hemocytes, a newly identified target tissue, involves binding of toxin to aminopeptidase N1 receptor](#). PLoS Pathog 17(3), e1009463, (2021), DOI: 10.1371/journal.ppat.1009463 (CellBrite® Red)

Metzger, W., et al. [Expansion and differentiation of human primary osteoblasts in two- and three-dimensional culture](#). Biotech Histochem 88(2), 86-102, (2013), DOI: 10.3109/10520295.2012.741262 (CellBrite® Blue)

Mohanty, S., et al. [Improving the efficacy of osteosarcoma therapy: combining drugs that turn cancer cell 'don't eat me' signals off and 'eat me' signals on](#). Mol Oncol 13(10), 2049-2061, (2019), DOI: 10.1002/1878-0261.12556 (CellBrite® Green)

Mohanty, S., et al. [Nanoparticle enhanced MRI can monitor macrophage response to CD47 mAb immunotherapy in osteosarcoma](#). Cell Death Dis 10(2), 36, (2019), DOI: 10.1038/s41419-018-1285-3 (CellBrite® Green)

Moquin-Beaudry, G., et al. [The Tumor-Immune Response Is Not Compromised by Mesenchymal Stromal Cells in Humanized Mice](#). J Immunol 203(10), 2735-2745, (2019), DOI: 10.4049/jimmunol.1900807 (CellBrite® NIR790)

- Morone, S., et al. [Overexpression of CD157 contributes to epithelial ovarian cancer progression by promoting mesenchymal differentiation](#). PLoS One 7(8), e43649, (2012), DOI: 10.1371/journal.pone.0043649 (CellBrite® Red)
- O'Neill, E., et al. [Imaging DNA Damage Repair In Vivo After \(177\)Lu-DOTATATE Therapy](#). J Nucl Med 61(5), 743-750, (2020), DOI: 10.2967/jnumed.119.232934 (CellBrite® Orange)
- O'Neill, J. D., et al. [Cross-circulation for extracorporeal support and recovery of the lung](#). Nature Biomedical Engineering 1(3), (2017), DOI: 10.1038/s41551-017-0037 (CellBrite® NIR790)
- Pajenda, G., et al. [Spatiotemporally limited BDNF and GDNF overexpression rescues motoneurons destined to die and induces elongative axon growth](#). Exp Neurol 261, 367-376, (2014), DOI: 10.1016/j.expneurol.2014.05.019 (CellBrite® Red)
- Prats-Mateu, B., et al. [Influence of HepG2 cell shape on nanoparticle uptake](#). Microsc Res Tech 77(8), 560-565, (2014), DOI: 10.1002/jemt.22374 (CellBrite® Orange)
- Shafa, M., et al. [Human induced pluripotent stem cell-derived lung progenitor and alveolar epithelial cells attenuate hyperoxia-induced lung injury](#). Cytotherapy 20(1), 108-125, (2018), DOI: 10.1016/j.jcyt.2017.09.003 (CellBrite® Orange, CellBrite® Red)
- Startek, J. B., et al. [Mouse TRPA1 function and membrane localization are modulated by direct interactions with cholesterol](#). Elife 8, (2019), DOI: 10.7554/eLife.46084 (CellBrite® Green)
- Suresh, M., et al. [In vitro model of inflammatory, hypoxia, and cancer stem cell signaling in pancreatic cancer using heterocellular 3-dimensional spheroids](#). bioRxiv, (2018), DOI: 10.1101/454397 (CellBrite® Green, CellBrite® Red, CellBrite® Blue, cell labeling before spheroid formation)
- Tarulli, E., et al. [Effectiveness of micron-sized superparamagnetic iron oxide particles as markers for detection of migration of bone marrow-derived mesenchymal stromal cells in a stroke model](#). J Magn Reson Imaging 37(6), 1409-1418, (2013), DOI: 10.1002/jmri.23897 (CellBrite® Green, CellBrite® Red)
- Tiptiri-Kourpeti, A., et al. [Lactobacillus casei Exerts Anti-Proliferative Effects Accompanied by Apoptotic Cell Death and Up-Regulation of TRAIL in Colon Carcinoma Cells](#). PLoS One 11(2), e0147960, (2016), DOI: 10.1371/journal.pone.0147960 (CellBrite® Red)
- Veraitch, O., et al. [Induction of hair follicle dermal papilla cell properties in human induced pluripotent stem cell-derived multipotent LNGFR\(+\)THY-1\(+\) mesenchymal cells](#). Sci Rep 7, 42777, (2017), DOI: 10.1038/srep42777 (CellBrite® Orange)
- Waehrens, L. N., et al. [Bovine lactadherin as a calcium-independent imaging agent of phosphatidylserine expressed on the surface of apoptotic HeLa cells](#). J Histochem Cytochem 57(10), 907-914, (2009), DOI: 10.1369/jhc.2009.953729 (CellBrite® Green)
- Weiss, T., et al. [NKG2D-Based CAR T Cells and Radiotherapy Exert Synergistic Efficacy in Glioblastoma](#). Cancer Res 78(4), 1031-1043, (2018), DOI: 10.1158/0008-5472.CAN-17-1788 (CellBrite® NIR790)
- Whitehead, J., et al. [Hydrogel mechanical properties are more important than osteoinductive growth factors for bone formation with MSC spheroids](#). bioRxiv, (2020), DOI: 10.1101/2020.03.10.986315 (CellBrite® NIR680)

- Winkel, B. M. F., et al. [Plasmodium sporozoites induce regulatory macrophages](#). PLoS Pathog 16(9), e1008799, (2020), DOI: 10.1371/journal.ppat.1008799 (CellBrite® Green)
- Yaekashiwa, N., et al. [Verification of Non-thermal Effects of 0.3–0.6 THz-Waves on Human Cultured Cells](#). Photonics 6(1), (2019), DOI: 10.3390/photonics6010033 (CellBrite® Green)
- Yamamoto, K., et al. [Development of a human neuromuscular tissue-on-a-chip model on a 24-well-plate-format compartmentalized microfluidic device](#). Lab Chip 21(10), 1897-1907, (2021), DOI: 10.1039/d1lc00048a (CellBrite® Green)
- Yamashita, A., et al. [Culture substrate-associated YAP inactivation underlies chondrogenic differentiation of human induced pluripotent stem cells](#). Stem Cells Transl Med 10(1), 115-127, (2021), DOI: 10.1002/sctm.20-0058 (CellBrite® Red)
- Yan, G., et al. [BET inhibition induces an anti-apoptotic adaptive response and therapeutic vulnerability to MCL1 inhibitors in breast cancer](#). bioRxiv, (2020), DOI: 10.1101/711895 (CellBrite® Orange)
- Yang, N. and Friedl, A. [Syndecan-1-Induced ECM Fiber Alignment Requires Integrin  \$\alpha\$ v \$\beta\$ 3 and Syndecan-1 Ectodomain and Heparan Sulfate Chains](#). PLoS One 11(2), e0150132, (2016), DOI: 10.1371/journal.pone.0150132 (CellBrite® Green)
- Yang, T., et al. [Exosome delivered anticancer drugs across the blood-brain barrier for brain cancer therapy in Danio rerio](#). Pharm Res 32(6), 2003-2014, (2015), DOI: 10.1007/s11095-014-1593-y (CellBrite® Red)
- Yu, M., et al. [In PC3 prostate cancer cells ephrin receptors crosstalk to beta1-integrins to strengthen adhesion to collagen type I](#). Sci Rep 5, 8206, (2015), DOI: 10.1038/srep08206 (CellBrite® Green)
- Zhang, Y., et al. [Evaluation of immunomodulatory effect of recombinant human granulocyte-macrophage colony-stimulating factor on polymorphonuclear cell from dogs with cancer in vitro](#). Vet Comp Oncol 15(3), 968-979, (2017), DOI: 10.1111/vco.12236 (CellBrite® Blue)
- Zhang, Y. and Yang, H. [A simple and robust fluorescent labeling method to quantify trophoblast fusion](#). Placenta 77, 16-18, (2019), DOI: 10.1016/j.placenta.2019.01.016 (CellBrite® Orange)
- Zulueta, A., et al. [Lung mesenchymal stem cells-derived extracellular vesicles attenuate the inflammatory profile of Cystic Fibrosis epithelial cells](#). Cell Signal 51, 110-118, (2018), DOI: 10.1016/j.cellsig.2018.07.015 (CellBrite® Green)
- Zulueta, A., et al. [Inflammatory role of extracellular sphingolipids in Cystic Fibrosis](#). Int J Biochem Cell Biol 116, 105622, (2019), DOI: 10.1016/j.biocel.2019.105622 (CellBrite® Green)

### **CellBrite® Fix Membrane Stains**

- Cross, K. L., et al. [Targeted isolation and cultivation of uncultivated bacteria by reverse genomics](#). Nat Biotechnol 37(11), 1314-1321, (2019), DOI: 10.1038/s41587-019-0260-6 (CellBrite® Fix 640 bacteria)
- Fabrice, T. N., et al. [Interactome and F-Actin Interaction Analysis of Dictyostelium discoideum Coronin A](#). Int J Mol Sci 21(4), (2020), DOI: 10.3390/ijms21041469 (CellBrite® Fix 640 bacteria, yeast particles)
- Ghilardi, S. J., et al. [Ventral Stress Fibers Induce Plasma Membrane Deformation in Human Fibroblasts](#). bioRxiv, (2021), DOI: 10.1101/2021.03.01.433420 (CellBrite® Fix 488)

Li, Q., et al. [Plasticity in the grazing ecophysiology of Florenciella \(Dichtyochophyceae\), a mixotrophic nanoflagellate that consumes Prochlorococcus and other bacteria](#). *Limnology and Oceanography* 66(1), 47-60, (2020), DOI: 10.1002/lno.11585 (CellBrite® Fix 488)

Malik, S. and Bahal, R. [Investigation of PLGA nanoparticles in conjunction with nuclear localization sequence for enhanced delivery of anti-miR phosphorothioates in cancer cells in vitro](#). *J Nanobiotechnology* 17(1), 57, (2019), DOI: 10.1186/s12951-019-0490-2 (CellBrite® Fix 488)

Mitchell, A., et al. [Human induced pluripotent stem cells ameliorate hyperoxia-induced lung injury in a mouse model](#). *Am J Transl Res* 12(1), 292-307, (2020), (CellBrite® Fix human iPSCs)

Santiana, M., et al. [Vesicle-Cloaked Virus Clusters Are Optimal Units for Inter-organismal Viral Transmission](#). *Cell Host Microbe* 24(2), 208-220 e208, (2018), DOI: 10.1016/j.chom.2018.07.006 (CellBrite® Fix 488, virus EV, SIM)

Switacz, V. K., et al. [Influence of Size and Cross-Linking Density of Microgels on Cellular Uptake and Uptake Kinetics](#). *Biomacromolecules* 21(11), 4532-4544, (2020), DOI: 10.1021/acs.biomac.0c00478 (CellBrite® Fix 640)

### **MemBrite® Fix Cell Surface Staining Kits**

De Ieso, M. L., et al. [Combined pharmacological administration of AQP1 ion channel blocker AqB011 and water channel blocker Bacopaside II amplifies inhibition of colon cancer cell migration](#). *Sci Rep* 9(1), 12635, (2019), DOI: 10.1038/s41598-019-49045-9 (MemBrite® Fix 488/513)

Delfarah, A., et al. [Inhibition of nucleotide synthesis promotes replicative senescence of human mammary epithelial cells](#). *J Biol Chem* 294(27), 10564-10578, (2019), DOI: 10.1074/jbc.RA118.005806 (MemBrite® Fix 640/660)

Egnuni, T., et al. [Evaluating the targeting efficiency of anti-EGFR functionalised nanoparticles to Head and Neck cancer cells for use in NIR-II optical window](#). *Preprints*, (2021), DOI: 10.20944/preprints202103.0687.v1 (MemBrite® Fix 594/615)

Graham, A. D., et al. [The development of a high throughput drug-responsive model of white adipose tissue comprising adipogenic 3T3-L1 cells in a 3D matrix](#). *Biofabrication* 12(1), 015018, (2019), DOI: 10.1088/1758-5090/ab56fe (MemBrite® Fix 640/660)

Le, P., et al. [Counting growth factors in single cells with infrared quantum dots to measure discrete stimulation distributions](#). *Nat Commun* 10(1), 909, (2019), DOI: 10.1038/s41467-019-08754-5 (MemBrite® Fix 640/660)

Mair, B., et al. [Scalable, FACS-Free Genome-Wide Phenotypic Screening](#). *bioRxiv*, (2019), DOI: 10.1101/612887 (MemBrite® Fix 488/515)

## CellBrite® and MemBrite® References for Specialized Applications

### Organoids, Spheroids, or Matrigel/Hydrogel Cultures

Anguiano, M., et al. [The use of mixed collagen-Matrigel matrices of increasing complexity recapitulates the biphasic role of cell adhesion in cancer cell migration: ECM sensing, remodeling and forces at the leading edge of cancer invasion](#). PLoS One 15(1), e0220019, (2020), DOI: 10.1371/journal.pone.0220019 (CellBrite® Red, cell labeling before Matrigel® culture)

Guimarães, C. F., et al. [3D flow-focusing microfluidic biofabrication: One-chip-fits-all hydrogel fiber architectures](#). Applied Materials Today 23, (2021), DOI: 10.1016/j.apmt.2021.101013 (CellBrite® Green, CellBrite Orange, cell labeling before hydrogel culture)

Suresh, M., et al. [In vitro model of inflammatory, hypoxia, and cancer stem cell signaling in pancreatic cancer using heterocellular 3-dimensional spheroids](#). bioRxiv, (2018), DOI: 10.1101/454397 (CellBrite® Green, CellBrite® Red, CellBrite® Blue, cell labeling before spheroid formation)

### Exosome and EV Labeling

Santiana, M., et al. [Vesicle-Cloaked Virus Clusters Are Optimal Units for Inter-organismal Viral Transmission](#). Cell Host Microbe 24(2), 208-220 e208, (2018), DOI: 10.1016/j.chom.2018.07.006 (CellBrite® Fix 488, virus EV, SIM)

### Super Resolution Imaging

Santiana, M., et al. [Vesicle-Cloaked Virus Clusters Are Optimal Units for Inter-organismal Viral Transmission](#). Cell Host Microbe 24(2), 208-220 e208, (2018), DOI: 10.1016/j.chom.2018.07.006 (CellBrite® Fix 488, virus EV, SIM)