CytoLiner™ Fixed Cell Membrane Stains

Novel lipophilic fluorescent dyes for selective staining of the plasma membrane in formaldehyde-fixed cells. Available in 6 colors, from blue to near-IR.



Product Description

CytoLinerTM Fixed Cell Membrane Dyes are novel lipophilic fluorescent dyes developed specifically for selective plasma membrane staining in fixed cells for microscopy. The dyes are engineered for robust and consistent staining of formaldehyde-fixed cells, and are suitable for downstream immunofluorescence staining protocols.

Novel Membrane Dyes Designed for Fixed Cells

While classic lipophilic membrane dyes like Dil can be used to stain formaldehyde-fixed cells, staining can be highly variable due to the poor solubility of the dyes. CytoLinerTM Dyes are a new generation of membrane dyes uniquely engineered to permit selective staining of the plasma membrane in fixed and mildly permeabilized cells. CytoLinerTM staining is compatible with formaldehyde fixation and mild detergent permeabilization before staining. Staining can tolerate blocking agents used during immunofluorescence staining protocols, although fish gelatin is recommended for best results, allowing subsequent staining with antibodies and other probes. The dyes also are compatible with poly-L-lysine coated cultureware and Transwell® permeable filter supports. CytoLinerTM Dyes are available in a selection of 6 colors from blue to near-IR.

Superior Staining Performance Over Classic Membrane Dyes Suitable for Co-Staining with Antibodies and Other Probes

Find the Right Stain for Your Application

CytoLinerTM Dyes are part of Biotium's family of unique and innovative cell membrane stains designed for specific applications. For workflows requiring staining live cells before fixation, we recommend our CellBrite® Fix Membrane Stains or MemBrite® Fix Cell Surface Staining Kits. Biotium also offers CF® Dye conjugated lectins, including Concanavalin A (Con A) and Wheat Germ Agglutinin (WGA), for visualizing cell surfaces in live or fixed cells. For longer term monitoring of cell surfaces in live cells, we recommend CellBrite® Steady Dyes. To find the right stain for your application, see our Membrane & Cell Surface Stains Comparison, or download our Membrane & Surface Stains Brochure.

Watch our video where Technical Applications Scientist II, Jacqueline Steenhuis PhD answers your top questions about Biotium's various membrane stains for fluorescence microscopy.

For additional support or product recommendations, contact us at techsupport@biotium.com.

CytoLiner™ Fixed Cell Membrane Stains

| Product | Ex/Em | Detection Channels | Size | Catalog Number |
|---------------------------------|------------|---------------------------|----------------|----------------|
| CytoLiner™ 410/450 | 406/446 nm | DAPI/Pacific Blue™ | 250 Labelings | <u>30131-T</u> |
| | | | 1000 Labelings | <u>30131</u> |
| CytoLiner™ 495/510 | 492/510 nm | FITC | 250 Labelings | <u>30132-T</u> |
| | | | 1000 Labelings | <u>30132</u> |
| CytoLiner™ 570/590 | 573/592 nm | Cy®3/TRITC | 250 Labelings | <u>30133-T</u> |
| | | | 1000 Labelings | <u>30133</u> |
| CytoLiner™ 650/675 | 647/674 nm | Cy®5 | 250 Labelings | <u>30134-T</u> |
| | | | 1000 Labelings | <u>30134</u> |
| CytoLiner™ 680/710 | 682/707 nm | Cy®5.5 | 250 Labelings | <u>30135-T</u> |
| | | | 1000 Labelings | <u>30135</u> |
| CytoLiner TM 785/815 | 787/819 nm | Alexa Fluor® 790 | 250 Labelings | <u>30140-T</u> |
| | | | 1000 Labelings | <u>30140</u> |

This datasheet was generated on November 3, 2025 at 02:14:54 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cytoliner-fixed-cell-membrane-stains/

Call us: 800-304-5357 Email: btinfo@biotium.com

Product attributes

| Probe cellular localization | Membrane/cell surface | | |
|-----------------------------|--|--|--|
| For live or fixed cells | For fixed cells | | |
| Assay type/options | Endpoint assay | | |
| Fixation options | Fix before staining (formaldehyde), Permeabilize before staining | | |
| Colors | Blue, Green, Red, Far-red, Near-infrared | | |