Oxazole Blue Homodimer (POPO™-1), 1 mM in DMSO

Oxazole Blue Homodimer, also known as POPOTM-1, is a blue-fluorescent, cell-impermeant nucleic acid stain that can be used to stain dead or fixed cells.

Product Description

Oxazole Blue Homodimer, also known as POPOTM-1, is a blue-fluorescent, cell-impermeant, high-affinity carbocyanine dimeric nucleic acid stain. The dye is essentially non-fluorescent in the absence of nucleic acids but exhibits strong excitation/emission maxima 433/457 nm when bound to nucleic acids. In addition, the dye is non-cytotoxic and may be used for long-term monitoring of viability in cell cultures.

- Dimeric nucleic acid stain
- Cell-impermeant
- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 433/457 nm
- Supplied at 1 mM in DMSO
- C41H54I4N6O2
- MW: 1170
- CAS number: 169454-15-3

See the table below for other chemical equivalents of Thermo Fisher Scientific's branded dead-cell selective nucleic acid dyes.

Biotium also offers unique NucSpot® Nuclear Stains for bright and specific nuclear staining in dead or fixed cells. The stains are available in a wide range of colors from green to near-IR. See our Cellular Stains Selection Guide and Cellular Stains Table for more information on other nuclear stains we offer.

| Product | Equivalent to | Color (Ex/Em) | Catalog No. |
|-----------------------------------------------|---------------|----------------------|--------------|
| Oxazole Blue, 1 mM in DMSO | PO-PROTM-1 | Blue (434/457 nm) | 40091 |
| Oxazole Blue Homodimer, 1 mM in DMSO | POPOTM-1 | Blue (433/457 nm) | 40093 |
| Oxazole Yellow, 1 mM in DMSO | YO-PRO®-1 | Green (491/506 nm) | 40089 |
| Oxazole Yellow Homodimer, 1 mM in DMSO | YOYO®-1 | Green (491/508 nm) | 40090 |
| TO lodide, 1 mM in DMSO | TO-PRO®-1 | Green (515/531 nm) | 40088 |
| Thiazole Orange Homodimer, 1 mM in DMSO | TOTO®-1 | Green (514/531 nm) | 40079 |
| Oxazole Red, 1 mM in DMSO | YO-PRO®-3 | Far-red (613/629 nm) | <u>40105</u> |
| Oxazole Red Homodimer, 1 mM in DMSO | YOYO®-3 | Far-red (612/631 nm) | 40106 |
| Thiazole Red, 1 mM in DMSO | TO-PRO®-3 | Far-red (642/657 nm) | 40087 |
| Thiazole Red Homodimer, 1 mM in DMSO | TOTO®-3 | Far-red (642/661 nm) | 40080 |

YOYO, YO-PRO, POPO, PO-PRO, TOTO, and TO-PRO are trademarks and registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on November 2, 2025 at 03:25:30 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/oxazole-blue-homodimer-popo-1-mm-in-dmso/

Product attribute

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

Apoptosis/viability marker
For live or fixed cells
Detection method/readout
Probe cellular localization
Assay type/options
Cell permeability
Colors
CAS number

Excitation/Emission