Oxazole Red (YO-PRO®-3), 1 mM in DMSO

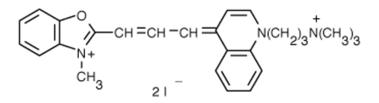
Bictum

Oxazole Red, also known as YO-PRO®-3, is a far-red fluorescent, cell-impermeant nucleic acid stain that can be used to stain dead or fixed cells.

Product Description

Oxazole Red, also known as YO-PRO®-3, is a far-red fluorescent, cell-impermeant nucleic acid stain that can be used to stain dead or fixed cells. The dye is essentially non-fluorescent in the absence of nucleic acids but exhibits strong excitation/emission maxima 613/629 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.

- Monomeric nucleic acid stain
- Cell-impermeant
- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 613/629 nm
- Supplied at 1 mM in DMSO
- C₂₆H₃₁I₂N₃O
- MW: 655



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

Biotium also offers unique <u>NucSpot® Nuclear Stains</u> 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 <u>Cellular Stains Selection Guide</u> and <u>Cellular Stains Table</u> for more information on other nuclear stains we offer.

Product	Equivalent to	Color (Ex/Em)	Catalog No.
Oxazole Blue, 1 mM in DMSO	PO-PRO™-1	Blue (434/457 nm)	<u>40091</u>
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)	<u>40089</u>
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)	<u>40088</u>
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)	<u>40087</u>
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.

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

Product attributes

Apoptosis/viability marker	Dead cell stain
For live or fixed cells	For fixed cells, For live/intact cells
Detection method/readout	Fluorescence microscopy, Flow cytometry
Probe cellular localization	Nucleus & cytoplasm
Assay type/options	No-wash staining, Real-time imaging
Cell permeability	Membrane impermeant
Colors	Far-red
CAS number	157199-62-7
Excitation/Emission	613/629 nm (with DNA)