

## Oxazole Red Homodimer (YOYO®-3), 1 mM in DMSO

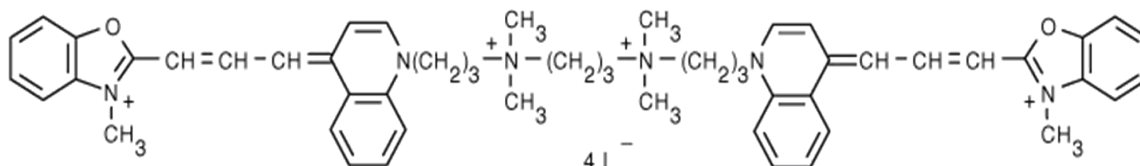
Oxazole Red Homodimer, also known as YOYO®-3, is a far-red fluorescent, cell-impermeant nucleic acid stain that can be used to stain dead or fixed cells.



### Product Description

Oxazole Red Homodimer, also known as YOYO®-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 612/631 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) = 612/631 nm
- Supplied at 1 mM in DMSO
- $C_{53}H_{62}N_4O_2$
- MW: 1323



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.
<a href="#">Oxazole Blue, 1 mM in DMSO</a>	PO-PRO™-1	Blue (434/457 nm)	<a href="#">40091</a>
<a href="#">Oxazole Blue Homodimer, 1 mM in DMSO</a>	POPO™-1	Blue (433/457 nm)	<a href="#">40093</a>
<a href="#">Oxazole Yellow, 1 mM in DMSO</a>	YO-PRO®-1	Green (491/506 nm)	<a href="#">40089</a>
<a href="#">Oxazole Yellow Homodimer, 1 mM in DMSO</a>	YOYO®-1	Green (491/508 nm)	<a href="#">40090</a>
<a href="#">TO Iodide, 1 mM in DMSO</a>	TO-PRO®-1	Green (515/531 nm)	<a href="#">40088</a>
<a href="#">Thiazole Orange Homodimer, 1 mM in DMSO</a>	TOTO®-1	Green (514/531 nm)	<a href="#">40079</a>
<a href="#">Oxazole Red, 1 mM in DMSO</a>	YO-PRO®-3	Far-red (613/629 nm)	<a href="#">40105</a>
<a href="#">Oxazole Red Homodimer, 1 mM in DMSO</a>	YOYO®-3	Far-red (612/631 nm)	<a href="#">40106</a>
<a href="#">Thiazole Red, 1 mM in DMSO</a>	TO-PRO®-3	Far-red (642/657 nm)	<a href="#">40087</a>
<a href="#">Thiazole Red Homodimer, 1 mM in DMSO</a>	TOTO®-3	Far-red (642/661 nm)	<a href="#">40080</a>

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

This datasheet was generated on June 15, 2026 at 06:22:58 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/oxazole-red-homodimer-yoyo-3-1-mm-in-dmso/>

### Product attributes

Apoptosis/viability marker	Dead cell stain
For live or fixed cells	For fixed cells, For live/f
Detection method/readout	Fluorescence microscop
Probe cellular localization	Nucleus & cytoplasm
Assay type/options	No-wash staining, Real
Cell permeability	Membrane impermeant
Colors	Far-red
CAS number	156312-20-8
Excitation/Emission	612/631 nm (with DNA)