Oxazole Yellow Homodimer (YOYO®-1), 1 mM in DMSO

nicture 4

Oxazole Yellow Homodimer, also known as YOYO®-1, is a green-fluorescent, cell-impermeant nucleic acid stain that can be used to stain dead or fixed cells.

Product Description

Oxazole Yellow Homodimer, also known as YOYO®-1, is a green-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 491/508 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) = 491/508 nm
- Supplied at 1 mM in DMSO
- C49H58I4N6O2
- MW: 1271

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 7, 2025 at 11:51:16 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/oxazole-yellow-homodimer-yoyo-1-1-mm-in-dmso/

Product attribute

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

Apoptosis/viability mark/
For live or fixed cells
Detection method/reado
Probe cellular localizatio
Assay type/options
Cell permeability
Colors
CAS number

Excitation/Emission