Glowing products for science

Oxazole Yellow (YO-PRO®-1), 1 mM in DMSO

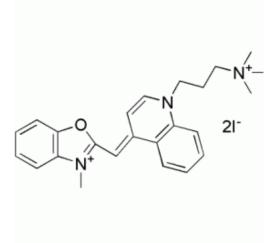
ulcitum

Oxazole Yellow, also known as YO-PRO®-1, is a green-fluorescent, cell-impermeant nucleic acid stain that can be used as a early marker of apoptosis.

Product Description

Oxazole Yellow, also known as YO-PRO®-1, is a green-fluorescent, cell-impermeant, high-affinity carbocyanine monomeric nucleic acid stain. It is essentially non-fluorescent in the absence of nucleic acids but exhibits excitation/emission maxima 491/506 nm when bound to nucleic acids. Oxazole Yellow is also used to identify apoptotic cells. Early apoptotic cells become permeant to Oxazole Yellow, but remain impermeant to propidium iodide (cat. no. 40016), a dead cell stain.

- Monomeric nucleic acid stain
- Cell-impermeant
- Early marker of apoptosis
- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 491/506 nm
- Supplied at 1 mM in DMSO
- C₂₄H₂₉I₂N₃O
- MW: 629



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.

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	Green		
CAS number	152068-09-2		
Excitation/Emission	491/506 nm (with DNA)		

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)	<u>40106</u>
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 09:20:48 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/oxazole-yellow-1mm-in-dmso/