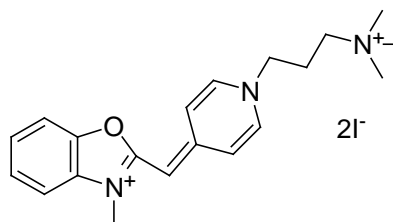


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Oxazole Blue, 1mM in DMSO**CATALOG NUMBER:** 40091**MOLECULAR INFORMATION:** C₂₀H₂₇I₂N₃O
Mwt: 579.26
CAS#: 157199-56-9**PROPERTIES:****Color & Form**

Yellow solution

Purity

≥ 95% by HPLC

 $\lambda_{ex}/\lambda_{em}$

~435/455nm (with DNA)

STORAGE AND HANDLING:

Store at -20°C and protect from light.

APPLICATION:

Oxazole blue, also known as PO-PRO-1, is a cell-impermeant, high-affinity carbocyanine monomeric nucleic acid stain that is essentially nonfluorescent in the absence of nucleic acids but exhibits excitation/emission maxima 435/455 nm when bound to nucleic acids. Oxazole blue is also used to identify apoptotic cells. Supplied at 1 mM in DMSO.

TOXICITY:

Unknown

FIRST AID:

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.