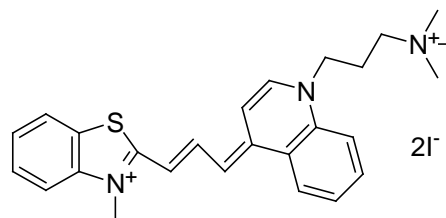


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Thiazole Red, 1mM in DMSO**CATALOG NUMBER:** 40087**MOLECULAR INFORMATION:** C₂₆H₃₁I₂N₃S
Mwt: 671.42**PROPERTIES:**

Color & Form Blue solution
Purity ≥ 95% by HPLC
λ_{ex}/λ_{em} ~642/661nm (with DNA)

STORAGE AND HANDLING:

Store at -20°C and protect from light.

APPLICATION:

Thiazole red, also known as TO-PRO-3, 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 ~642/661 nm when bound to nucleic acids. Thiazole red is also used as a nuclear counterstain and dead cell indicator.

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.