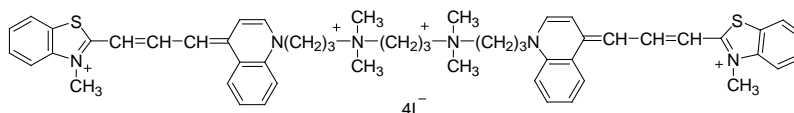


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Thiazole Red Homodimer, 1 mM in DMSO**CATALOG #** 40080**MOLECULAR INFORMATION:** C₅₃H₆₂I₄N₆S₂
MWt: 1354.85**PROPERTIES:****Color & Form** Blue solution
Absorption/Emission 642/660 nm (DNA-bound)**STORAGE AND HANDLING:** Stored desiccated at 4 °C and protect from light**APPLICATION:** Thiazole red homodimer (TRhD), also known as TOTO-3, is a cell-impermeant, high-affinity nucleic acid stain based upon a symmetric cyanine dye dimer that is essentially nonfluorescent in the absence of nucleic acids that exhibits excitation/emission maxima ~642/660 nm when bound to nucleic acids.**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.