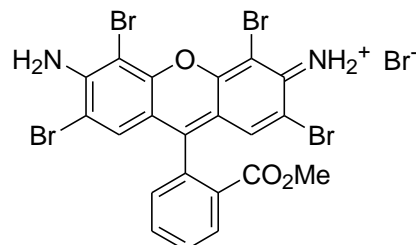


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 2', 4', 5' 7' - tetrabromorhodamine 123 bromide

CATALOG NUMBER: 70013

MOLECULAR INFORMATION: C₂₁H₁₃Br₅N₂O₃
Mwt: 740.86



PROPERTIES:

Color & Form	Orange solid
Purity	≥ 95% by HPLC
Solubility	Soluble in DMSO
$\lambda_{\text{ex}} \backslash \lambda_{\text{em}}$	524nm \ 550nm (MeOH)
Extinction Coefficient (ϵ)	>85,000

STORAGE AND HANDLING:

Store at 4 °C and protect from light

APPLICATION:

Tetrabromorhodamine 123 can be used to generate singlet oxygen (¹O₂) in mitochondrial membranes^{1,2} The dye has a quantum yield for singlet oxygen generation of 0.65-0.7 and is particularly toxic to carcinoma cells.³

Ref.: 1) *Photochem. Photobiol.* **55**, 81(1992); 2) *Proc. SPIE* **997**, 48 (1988); 3) *Cancer Res.* **49**, 3961(1989).

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.