

Tetrabromorhodamine 123, bromide

Tetrabromorhodamine 123 can be used to generate singlet oxygen (1O_2) 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.



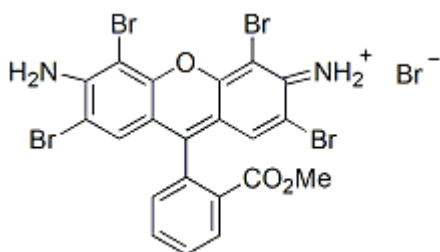
Product attributes

Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining, Real-time imaging
Colors	Red
Excitation/Emission	524/550 nm

Product Description

Tetrabromorhodamine 123 can be used to generate singlet oxygen (1O_2) 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

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 524/550 \text{ nm}$
- $\epsilon (\text{MeOH}) = 91,000$
- Orange red solid soluble in DMSO or DMF
- Store at 4 °C and protect from light, especially in solution
- $C_{21}H_{13}Br_5N_2O_3$
- MW: 740.86



References

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