Rhodamine 123

Rhodamine 123 is a popular green fluorescent mitochondrial dye that stains mitochondria in living cells in a membrane potential-dependent fashion. It is widely used in flow cytometry



Product Description

Rhodamine 123 is a popular green fluorescent mitochondrial dye that stains mitochondria in living cells in a membrane potential-dependent fashion. It is widely used in flow cytometry studies involving mitochondrial membrane potential. Please also see Dihydrorhodamine 123 and its HCl salt (10055 and 10056).

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 505/534 nm
- ϵ (505 nm, MeOH) = 97,000
- Orange red solid soluble in DMSO or DMF
- Store at 4°C and protect from light, especially in solution
- C₂₁H₁₇CIN₂O₃
- MW: 381
- [62669-70-9]

References

- 1. Cytometry 17, 50 (1994).
- 2. Science 218, 1117 (1982).
- 3. J Cell Biol 88, 526 (1981).

Assay type/options Real-time imaging studies involving mitochondrial membrane potential. Cell permeability Membrane permeant Apoptosis/viability marker Mitochondrial potential Potential dependence Mitochondrial potential-dependent

> Colors Red Excitation/Emission 505/534 nm

Product attributes

Probe cellular localization

For live or fixed cells

CAS number

Call us: 800-304-5357 Email: btinfo@biotium.com

62669-70-9

Mitochondria

For live/intact cells

This datasheet was generated on November 11, 2025 at 07:29:20 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rhodamine-123/