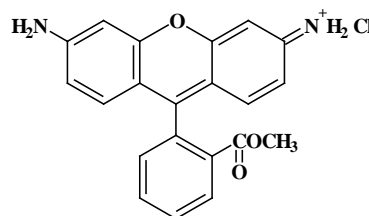


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Rhodamine 123**CATALOG NUMBER:** 70010**MOLECULAR INFORMATION:** C₂₁H₁₇ClN₂O₃
Mwt: 381
[62669-70-9]**PROPERTIES:**

| | |
|---|---|
| Color & Form | Orange red solid |
| Purity | ≥ 96% by HPLC |
| Solubility | Soluble in DMSO and DMF |
| $\lambda_{\text{ex}} \backslash \lambda_{\text{em}}$ | 505nm \ 534nm (in MeOH) |
| Extinction Coefficient (ϵ) | 97,000 M ⁻¹ cm ⁻¹ (505 nm; in MeOH) |

STORAGE AND HANDLING:

Stored desiccated at 4 °C upon receipt. Protect from light, especially when in solution.

APPLICATION:

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

Ref: 1) Millot, J.M., et al. *Cytometry* **17**, 50(1994); 2) Bernal, S.D., et al *Science* **218**, 1117(1982); 3) Johnson, L.V., et al. *J. Cell Biol.* **88**, 526(1981).**TOXICITY:**

Reproductive Effects: TDLo: 60 mg/kg intraperitoneal, mouse (7-10D preg) (post implantation mortality)

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.