

## Product Information

### Rhodamine 123

Cat. No.	Product Name	Unit Size
C044	Rhodamine 123	25 mg

**Storage upon receipt:**

- -20°C
- Protect from light

#### Product Description

**Rhodamine 123** is a cell-permeant, cationic, green-fluorescent dye that is readily sequestered by active mitochondria without cytotoxic effects. This product has been used to assay mitochondrial membrane potential in populations of apoptotic cells..

#### Specifications:

**Chemical name:** Rhodamine 123

**Molecular Formula:** C<sub>21</sub>H<sub>17</sub>ClN<sub>2</sub>O<sub>3</sub>

**Molecular Weight:** 380.83

**CAS Number:** 62669-70-9