## Dihydrorhodamine 123, Dihydrochloride Salt

atóstum

Dihydrorhodamine 123 dihydrochloride is functionally equivalent to dihydrorhodamine 123 (10055) but with increased stability toward air oxidation and light during storage.

## **Product Description**

Dihydrorhodamine 123 dihydrochloride is functionally equivalent to dihydrorhodamine 123 (10055) but with increased stability toward air oxidation and light during storage.

- White solid soluble in DMF
- Store at -20°C and protect from light, especially in solution
- C<sub>20</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>3</sub>
- MW: 419

2HCI

## References

- 1. Br J of Pharmacol (2010) doi: 10.1111/j.1476-5381.2010.01120.x
- 2. J Immunol Meth178, 89 (1995).
- 3. Biochemistry 34, 3544 (1995).
- 4. Eur. J. Biochem. 217, 973 (1993).
- 5. Arc. Biochem. Biophys. 302, 348 (1993).
- 6. Biometals (2015), doi: 10.1007/s10534-015-9873-5

This datasheet was generated on November 7, 2025 at 03:25:47 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dihydrorhodamine-123-dihydrochloride-salt/

## **Product attributes**

Call us: 800-304-5357

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	505/534 nm (end product)

Email: btinfo@biotium.com