

Dihydrorhodamine 123, Dihydrochloride Salt

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



Product attributes

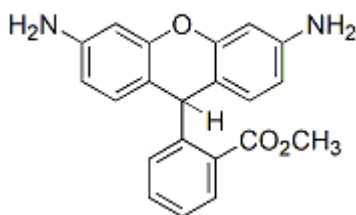
Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining, Real-time imaging
Colors	Red
Excitation/Emission	505/534 nm (end product)

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
- C₂₀H₂₀Cl₂N₂O₃
- MW: 419

2HCl



References

1. Br J of Pharmacol (2010) doi: 10.1111/j.1476-5381.2010.01120.x
2. J Immunol Meth 178, 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](https://doi.org/10.1007/s10534-015-9873-5)

This datasheet was generated on June 11, 2026 at 12:54:28 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/dihydrorhodamine-123-dihydrochloride-salt/>