

Dihydroethidium (Hydroethidium)

Dihydroethidium (also called hydroethidium) is the chemically reduced form of the commonly used DNA dye ethidium bromide. The probe is useful to detect oxidative activities in viable cells, including respiratory burst in phagocytes.



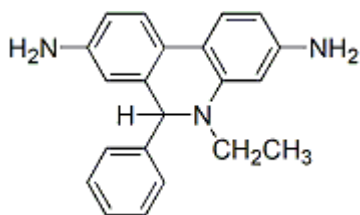
Product attributes

CAS number	38483-26-0
Excitation/Emission	518/605 nm (end product)

Product Description

Dihydroethidium (also called hydroethidium) is the chemically reduced form of the commonly used DNA dye ethidium bromide. The probe is useful to detect oxidative activities in viable cells, including respiratory burst in phagocytes. Dihydroethidium itself has blue fluorescence ($\lambda_{Ex}/\lambda_{Em}$: 355/420 nm) in cells, while the oxidized form ethidium has red fluorescence ($\lambda_{Ex}/\lambda_{Em}$: 518/605 nm) upon DNA intercalation.

- Off-white to light brown solid soluble in DMF or DMSO
- Store at -20°C under nitrogen or argon, and protect from light, especially when in solution
- C₂₁H₂₁N₃
- MW: 315
- [38483-26-0]



References

1. J Applied Technology (2010) DOI 10.1002/jat.1599
2. J Immunol Meth 170, 117 (1994).
3. FEMS Microbiol Lett 122, 187 (1994).
4. FEMS Microbiol Lett 101, 173 (1992).
5. J Histochem Cytochem 34, 1109 (1986).

This datasheet was generated on June 6, 2026 at 08:52:11 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dihydroethidium-also-called-hydroethidium/>