DiSC2(5) (High Purity)

DiSC₂(5) (3,3'-Diethylthiadicarbocyanine, iodide) binds to DNA-DNA or PNA-DNA duplexes with a large blue-shift of the absorption spectrum. Binding of the dye to PNA-DNA hybrid, in particular, blue-shifts the dye absorption spectrum by as much as 115 nm, offering a visual method to detect nucleic acid hybridization (1,2).

sidum 80

Product attributes

CAS number

514-73-8

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

DiSC₂(5) (3,3'-Diethylthiadicarbocyanine, iodide) binds to DNA-DNA or PNA-DNA duplexes with a large blue-shift of the absorption spectrum. Binding of the dye to PNA-DNA hybrid, in particular, blue-shifts the dye absorption spectrum by as much as 115 nm, offering a visual method to detect nucleic acid hybridization (1,2).

- λ_{Abs} (pH 7.5 buffer) = 646 nm (free dye); ~586 nm (dsDNA); ~531 nm (PNA-DNA)
- Dark solid soluble in DMSO or MeOH
- Store at 4°C
- C₂₃H₂₃IN₂S₂
- MW: 518.48
- [514-73-8]

References

- 1. Nucleic Acid Res 30(2), 3 (2002).
- 2. J Am Chem Soc 121, 2686 (1999).

This datasheet was generated on November 4, 2025 at 10:31:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/disc25-33-diethylthiadicarbocyanine-iodidehigh-purity/