

DODC (High Purity)

DODC (DiOC₂(5)) has been reported to bind to triplex DNA as well as dimeric hairpin quadruplex DNA (1). Binding of the dye to triplex DNA quenches the fluorescence of the dye and red-shifts the dye absorption peak.

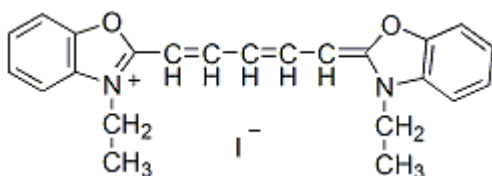


Product attributes

Product Description

DODC (DiOC₂(5)) has been reported to bind to triplex DNA as well as dimeric hairpin quadruplex DNA (1). Binding of the dye to triplex DNA quenches the fluorescence of the dye and red-shifts the dye absorption peak.

- $\lambda_{Abs} = 582 \text{ nm}$
- Red solid soluble in DMSO or MeOH
- Store at 4 °C
- C₂₃H₂₃N₂O₂
- MW: 486.35



References

1. J Am Chem Soc 122, 424 (2000).
2. PNAS 93, 2635 (1996).

This datasheet was generated on June 23, 2026 at 10:00:25 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/dodc-dioc25high-purity/>