

Product Information

DODC, (DiOC₂(5)) (high purity)

Catalog Number: 40040

Unit Size: 50mg

Storage and Handling:

Store at 4°C. Guaranteed for up to one year from the date of receipt when stored properly.

Spectral Properties:

λ_{abs} = 582nm

Product Description:

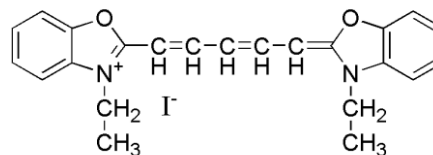
Molecular Formula: C₂₃H₂₃IN₂O₂

Molecular Weight: 486.35

Color and Form: Red solid

Solubility: Soluble in DMSO or MeOH

Application: DODC has been reported to bind to triplex DNA as well as dimeric hairpin quadruplex DNA.^{1,2} Binding of the dye to triplex DNA quenches the fluorescence of the dye and red-shifts the dye absorption peak.



References

- 1) *J. Am. Chem. Soc.* **122**, 424 (2000)
- 2) *Proc. Natl. Acad. Sci. USA* **93**, 2635 (1996)