

## DODC (High Purity)

DODC (DiOC<sub>2</sub>(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.

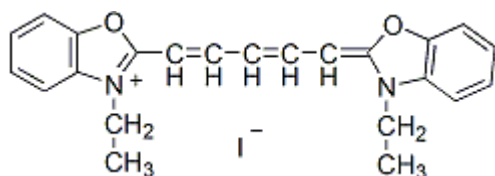


### Product attributes

## Product Description

DODC (DiOC<sub>2</sub>(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
- C<sub>23</sub>H<sub>23</sub>IN<sub>2</sub>O<sub>2</sub>
- MW: 486.35



## References

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

This datasheet was generated on December 30, 2025 at 05:56:55 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dodc-dioc25high-purity/>