

DiOC6(3)

DiOC₆(3) (3,3'-Dihexyloxacarboyanine,iodide) is a green fluorescent membrane dye that has been used to stain the ER in both live and fixed cells. The dye has also been used to study structural interactions and dynamics of the ER in neurons and yeast.



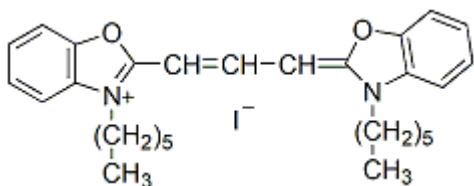
Product attributes

CAS number	53213-82-4
Probe cellular localization	Endoplasmic reticulum
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Potential dependence	Slow response (translational) membrane potential dye
Colors	Green
Excitation/Emission	484/501 nm

Product Description

DiOC₆(3) (3,3'-Dihexyloxacarboyanine,iodide) is a green fluorescent membrane dye that has been used to stain the ER in both live and fixed cells (1,2). The dye has also been used to study structural interactions and dynamics of the ER in neurons (3,4) and yeast (5). In addition, the dye is widely used as a membrane potential dye (6).

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 484/501 nm
- Orange solid soluble in DMSO or DMF
- Store at 4 °C and protect from light, especially in solution
- C₂₉H₃₇N₂O₂
- MW: 572.53
- [53213-82-4]



References

1. J Cell Biol 103, 1557 (1986).
2. Cell 38, 101 (1984).
3. J Cell Biol 127, 1021 (1994).
4. Nature, 310, 53 (1984).
5. Cell Motil Cytoskel 25, 111 (1993).

This datasheet was generated on June 15, 2026 at 01:29:35 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dioc63-33-dihexyloxacarboyanineiodide/>