

DiO

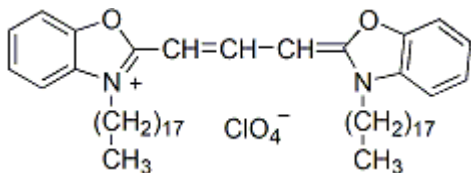
DiO, also called DiOC₁₈(3), is a green fluorescent, lipophilic carbocyanine dye that is widely used as a lipophilic tracer.



Product Description

DiO, also called DiOC₁₈(3), is a green fluorescent, lipophilic carbocyanine dye that is widely used as a lipophilic tracer. This membrane dye has applications similar to those of Dil (#60010). However, DiO has green fluorescent emission and the lateral diffusion rate on the membranes is generally slower than that of Dil. DiO and Dil are often used together in dual color studies.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 484/501 nm
- ϵ (MeOH) = 150,000
- Yellow orange solid soluble (with heating) at 2 mM in DMF, 1 mM in DMSO, or 2 mM in 1:1 DMSO:ethanol
- Store at 4 °C and protect from light, especially in solution
- C₅₃H₈₅ClN₂O₆
- MW: 882
- Cas number: 34215-57-1
- Cas name: Benzoxazolium, 3-octadecyl-2-[3-(3-octadecyl-2(3H)-benzoxazolylidene)-1-propenyl]-, perchlorate



Note: DiO has low solubility, a tendency to form aggregates and slow lateral diffusion rate. See our DiO alternative [Neuro-DiO \(60015\)](#), which is more soluble in membranes and does not form non-fluorescent aggregates. We also offer [CellBrite™ Green \(30021\)](#), a ready-to-use Neuro-DiO solution for cell membrane labeling, and [Neuro-DiO in Vegetable Oil \(60019\)](#) for microinjection studies.

References

1. Biochemistry 19, 6050 (1980)
2. J. Immunology 127, 893 (1981)
3. Trends Neurosci 9, 333 (1989)
4. J Cell Biol 103, 171 (1986)

This datasheet was generated on June 14, 2026 at 07:05:46 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dio-33-dioctadecyloxacarbocyanine-perchlorate/>

Product attributes

CAS number	34215-57-1
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	Co-cultures, Extended staining (several days to weeks)
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
Colors	Green
Excitation/Emission	484/501 nm