

DiOC16(3)

DiOC16(3) is a green fluorescent lipophilic cyanine dye with two C16 hydrocarbon chains that can be used to stain cytoplasmic membranes.



Product attributes

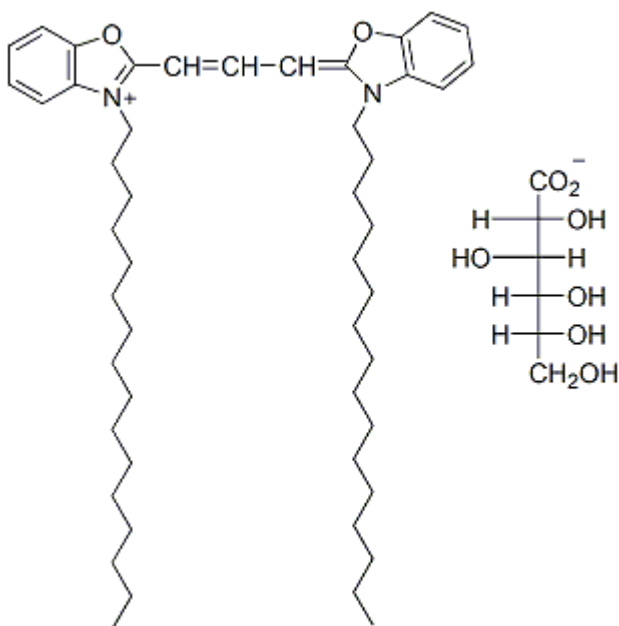
CAS number	161433-32-1
For live or fixed cells	For live/intact cells
Colors	Green
Excitation/Emission	484/501 nm

Product Description

DiOC16(3) is a green fluorescent lipophilic cyanine dye with two C16 hydrocarbon chains. The dye can be used to stain cytoplasmic membranes.

DiOC16 has been combined with DPA (cat # 60037) to detect cell membrane potential change with high sensitivity. See the [DiO/DPA Membrane Potential Detection Kit \(30037\)](#) for more information).

- Orange solid soluble in DMF and DMSO at >5 mg/mL
- $\lambda_{Abs}/\lambda_{Em}$ (MeOH) = 484/501 nm
- ϵ = ~130,000
- Store at 4 °C and protect from light
- $C_{55}H_{88}N_2O_9$
- MWt: 921.30
- [161433-32-1]



This datasheet was generated on June 5, 2026 at 12:13:13 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dioc163/>