

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
- $\lambda_{\text{Abs}}/\lambda_{\text{Em}}$ (MeOH) = 484/501 nm
- ϵ = ~130,000
- Store at 4 °C and protect from light
- $\text{C}_{55}\text{H}_{88}\text{N}_2\text{O}_9$
- MWt: 921.30
- [161433-32-1]

