

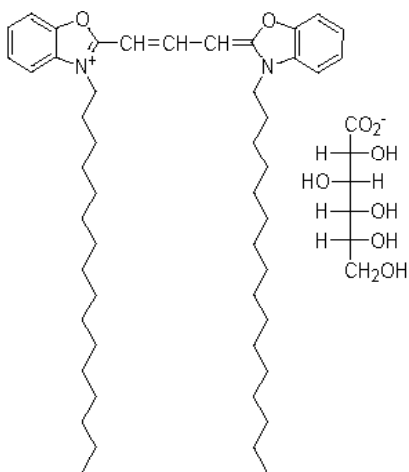
Product Information

DiOC16(3)

Catalog Number: 60038

Packaging Size: 25 mg

Molecular Information:

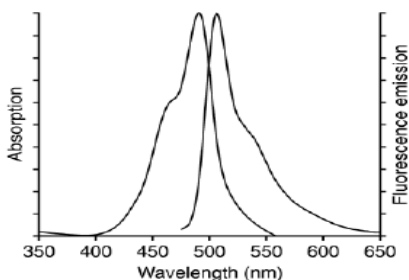


Mol. formula: C₅₅H₈₈N₂O₉; Mwt: 921.30

Color and Form: Orange solid

Solubility: Soluble in DMF or DMSO at >5 mg/mL (Note: sonication or heating may be applied to facilitate dissolution). Solution can be stored at room temperature, but must be protected from light.

Spectral Properties: $\lambda_{abs}/\lambda_{em}$ (MeOH) = 484/501 nm; ϵ = ~130,000



Storage and Handling

Store at room temperature or below. Shelf-life is at least 1 year from the date of receipt.

Product Description

DiOC16(3) (full chemical name: 3,3'-dihexadecyloxycarbocyanine) is a lipophilic cyanine dye with two C16 hydrocarbon chains. It can be optimally excited by the 488 nm argon laser line with green fluorescence emission (ex/em = 484/501 nm). The dye can be used to stain cytoplasmic membranes. Recently, DiOC16 has been combined with DPA (cat # 60037) to detect cell membrane potential change with high sensitivity (Bradley, et al. J. Neurosci. 29(29), 9197(2009)). Please also see our DiO/DPA membrane potential detection kit (cat# 30037).

Reference:

- Bradley, et al. J. Neurosci. 29(29), 9197(2009).

Safety

See MSDS.

First aid

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Other related products

- DiO/DPA membrane potential detection kit (cat# 30037).
- Dipicrylamine (Cat# 60037).
- Di-4-ANEPPS (Cat# 61010)
- SynpatoGreen C4 (or FM1-43) (Cat# 70020)
- Tetrodotoxin (Cat# 00061, 00062)
- Fluo-3, AM (Cat# 50014)

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.