

Product Information

DiOC₆(3)

Catalog Number	Unit Size
C050	25 mg

Storage upon receipt:

- -20°C
- Protect from light

Product Description

DiOC₆(3) is a cell-permeant, green-fluorescent, lipophilic dye that is selective for the mitochondria of live cells, when used at low concentrations. At higher concentrations, the dye may be used to stain other internal membranes, such as the endoplasmic reticulum.

Specifications:

Chemical Name: 3,3'-dihexyloxacarbocyanine iodide

Molecular Formula: C₂₉H₃₇IN₂O₂

Molecular Weight: 572.53

CAS Number: 53213-82-4