

DiOC6, Intracellular probe ab189808

画像数 1

製品の概要

製品名	DiOC6, Intracellular probe
製品の概要	Cell-permeant lipophilic dye for staining internal membranes (e.g., mitochondrial, Golgi membranes) of live & fixed cells. Concentration-dependent staining.
特記事項	<p>Biological Description: Cell-permeant, lipophilic dye that stains internal membranes of both live and fixed cells, such as the mitochondrial, Golgi and endoplasmic reticulum membranes. Staining patterns are concentration dependent.</p> <p>Purity: &gt;95%</p> <p>Chemical name: 3,3'-Dihexyloxacarbocyanine Iodide</p> <p>[CAS]:53213-82-4</p> <p>Molecular Formula:C<sub>29</sub>H<sub>37</sub>IN<sub>2</sub>O<sub>2</sub></p> <p>Molecular Weight: 572.53</p> <p>Excitation: 484</p> <p>Emission: 501</p> <p>Storage: room temperature, desiccated, protect from light</p> <p>Solubility: DMSO</p>
アプリケーション	適用あり: Other

製品の特性

保存方法Store at +4°C. Please refer to protocols.

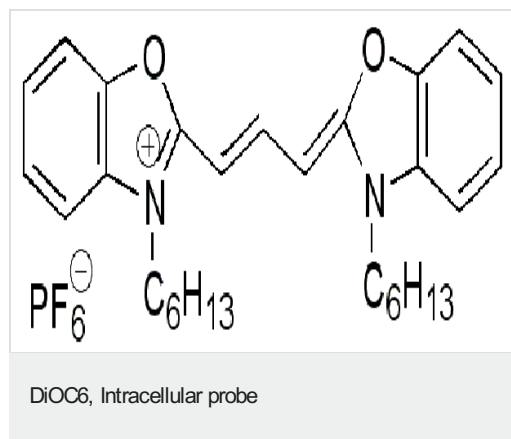
内容	100 mg
DiOC6 Intracellular probe	1 x 100mg

アプリケーション

The Abpromise guarantee      Abpromise保証は、次のテスト済みアプリケーションにおけるab189808の使用に適用されます  
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
Other		Use at an assay dependent concentration. Stains internal membranes of Golgi, ER & Mitochondria - Conc. dependent

## 画像



**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

## Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors