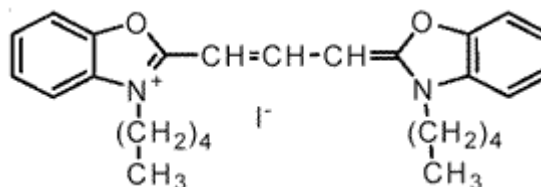


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** DiOC<sub>5</sub>(3), Iodide (3,3'-Diethyloxycarbocyanine, iodide)**CATALOG #:** 70007**MOLECULAR INFORMATION:** C<sub>27</sub>H<sub>33</sub>IN<sub>2</sub>O<sub>2</sub>  
Mwt: 544.47**PROPERTIES:**

**Color & Form** Orange solid  
**Purity** ≥ 98% TLC  
**Solubility** Soluble in DMF or DMSO  
**Absorption/Emission** 482/497 nm (in MeOH)  
**Extinction Coefficient** 120,000 cm<sup>-1</sup> M<sup>-1</sup> (in MeOH)

**STORAGE AND HANDLING:**

Store at 4°C and protect from light, especially when in solution.

**APPLICATION:**DiOC<sub>5</sub>(3) and DiOC<sub>6</sub>(3) are two of the most widely used carbocyanine dyes used for membrane potential measurements.

Ref: 1) J. Biol. Chem. 270, 3788 (1995).

**TOXICITY:**

Unknown

**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.

**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.