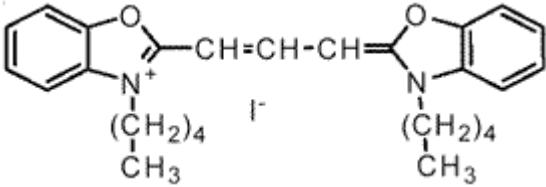


3,3'-Dipentylloxacarboyanine, iodide

Instruction Manual

| | |
|---------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog Number | PK-CA707-70007 |
| Description | DiOC ₅ (3) and DiOC ₆ (3) are two of the most widely used carbocyanine dyes used for membrane potential measurements. |
| Quantity | 100 mg |
| Excitation / Emission Maxima | $\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 482/497 nm Extinction Coefficient: 120,000 cm ⁻¹ M ⁻¹ (in MeOH) |
| Molecular Structure |  |
| Molecular Weight / Molecular Formula | 544.47 Da; C ₂₇ H ₃₃ IN ₂ O ₂ |
| Purity | ≥98% (TLC) |
| Appearance / Formulation / Solubility | Orange solid; soluble in DMSO or DMF. |
| Storage & Stability | Store at -20°C in a dry environment. Protect from light. |
| Applications | See Description |
| References | J. Biol. Chem. 270, 3788 (1995) |
| Caution | 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. |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.