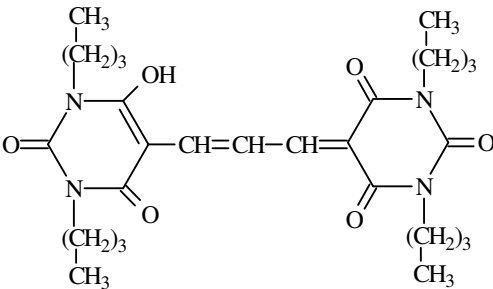


Instruction Manual

Catalog Number	PK-CA707-61011
Description	DiBAC ₄ (3) is a Bis-oxanol type so-called translational membrane potential dye that redistributes within the cell membrane when membrane potential changes. The fluorescence intensity of the dye is enhanced when the dye enters the cell membrane as a result of membrane depolarization. The rate of fluorescence response of the dye to membrane potential change is usually slower than that of the styryl dye di-4-ANEPPS, but the fluorescence signal change for the former is significantly larger than for the latter dye. Since an argon laser (488 nm) can be used for DiBAC ₄ (3) excitation, it is applicable for flow cytometry and confocal microscopy. A new and potentially very important application for DiBAC ₄ (3) is its use in high throughput drug screening.
Quantity	25 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 493/516 nm; Extinction coefficient = 123,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	516 Da; C ₂₇ H ₄₀ N ₄ O ₆
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Orange solid; soluble in DMSO and EtOH.
Storage & Stability	Store desiccated at $\leq 4^{\circ}\text{C}$. Protect from light, especially when in solution.
Applications	See Description
References	<ol style="list-style-type: none"> 1) Chem. Phys. Lipids 69, 137(1994) 2) Biochim. Biophys. Acta 771, 208(1984) 3) J. Biomol. Screening 1(2), 75(1996)
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.