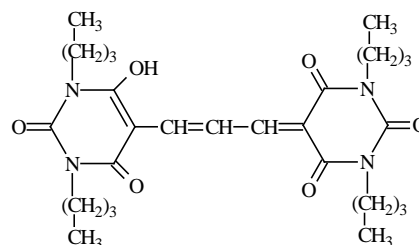


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** DiBAC<sub>4</sub>(3) ( bis-(1,3-dibarbituric acid)-trimethine oxanol)**CATALOG #:** 61011**MOLECULAR INFORMATION:** C<sub>27</sub>H<sub>40</sub>N<sub>4</sub>O<sub>6</sub>  
Mwt: 516  
[70363-83-6]**PROPERTIES:**

<b>Color &amp; Form</b>	Orange solid.
<b>Purity</b>	≥ 99% by HPLC
<b>Solubility</b>	Soluble in DMSO or EtOH
$\lambda_{ex} / \lambda_{em}$	493 nm / 516 nm (MeOH)
<b>Extinction Coefficient</b>	123,000 (MeOH)

**STORAGE AND HANDLING:**

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

**APPLICATION:**

DiBAC<sub>4</sub>(3) is a so-called translational membrane potential dye that redistributes within the cell membrane when membrane potential changes.<sup>1,2</sup> The fluorescence 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. A new and potentially very important application for DiBAC<sub>4</sub>(3) is its use in high throughput drug screening.<sup>3</sup>

Ref: 1) *Chem. Phys. Lipids* **69**, 137(1994); 2) *Biochim. Biophys. Acta* **771**, 208(1984); 3) *J. Biomol. Screening* **1**(2), 75(1996).**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.