DiBAC4(3)

 $\mathsf{DiBAC4}(3)$ is a translational membrane potential dye that redistributes within the cell membrane when membrane potential changes.



Product attributes

CAS number	70363-83-6
For live or fixed cells	For live/intact cells
Potential dependence	Slow response (translational) membrane potential dye
Colors	Green
Excitation/Emission	403/516 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

DiBAC4(3) (Bis-(1,3-Dibarbituric acid)-trimethine oxanol) is a so-called translational membrane potential dye that redistributes within the cell membrane when membrane potential changes (1,2). 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. DiBAC4(3) has applications in high-throughput drug screening.

- λεx/λεm (MeOH) = 493/516 nm
- Orange solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- C₂₇H₄₀N₄O₆
- MW: 519
- [70363-83-6]

References

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

This datasheet was generated on November 15, 2025 at 05:26:52 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dibac43-bis-13-dibarbituric-acid-trimethine-oxanol/