

DiBAC4(3)

DiBAC₄(3) is a translational membrane potential dye that redistributes within the cell membrane when membrane potential changes.



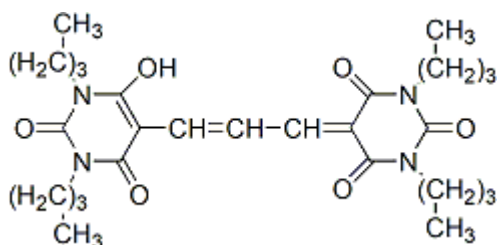
Product attributes

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	493/516 nm

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.

- $\lambda_{Ex}/\lambda_{Em}$ (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 June 19, 2026 at 09:55:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dibac43-bis-13-dibarbituric-acid-trimethine-oxanol/>