Di-12-ANEPPQ

Di-12-ANEPPQ is a fast-responding membrane potential dye. It is more hydrophobic than Di-8-ANEPPQ and is useful for potential-sensitive retrograde labeling of neurons.



Product attributes

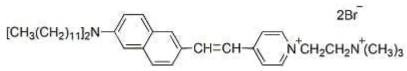
For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	522/720 nm (see product description)

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

Product Description

Di-12-ANEPPQ is a fast-responding membrane potential dye. It is more hydrophobic than Di-8-ANEPPQ and is useful for potential-sensitive retrograde labeling of neurons.

- λεx/λεm (MeOH) = 522/720 nm*
- Orange red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light, especially in solution
- C46H75Br2N3
- MW: 827.4



* Note that the spectra of styryl dyes generally undergo a large blue shift in lipids compared to organic solvent. See the spectra for styryl dyes in liposomes under product images.

References

- 1. J Neurosci Meth 70, 121 (1996).
- 2. J Neurosci Meth 70, 111 (1996).

This datasheet was generated on November 3, 2025 at 08:58:13 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/di-12-aneppg/