

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.



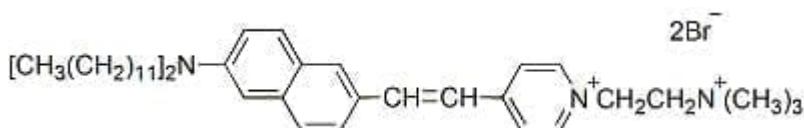
Product attributes

| | |
|-------------------------|--------------------------------------|
| 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) |

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.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 522/720 nm*
- Orange red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light, especially in solution
- $C_{46}H_{75}Br_2N_3$
- 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 January 2, 2026 at 05:38:37 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/di-12-aneppq/>