

Di-4-ANEPPS

Di-4-ANEPPS is a fast-responding membrane potential dye. Changes in the membrane potential of the cell correlate with changes in the fluorescence excitation intensity of the dye.



Product attributes

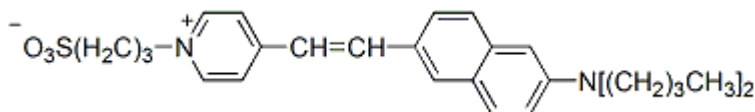
CAS number	90134-00-2
For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	496/705 nm (see product description)

Product Description

Di-4-ANEPPS is a fast-responding membrane potential dye (1). Changes in the membrane potential of the cell correlate with changes in the fluorescence excitation intensity of the dye. Di-4-ANEPPS is internalized in cells rapidly, so it is mainly used for short-term experiments. We also offer [Di-8-ANEPPS](#) (catalog number [61012](#)), which is more lipophilic and better retained in the outer leaflet of the plasma membrane.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 496/705 nm*
- Orange red solid soluble in DMF, ethanol or DMSO
- Store at 4 °C and protect from light, especially in solution
- $C_{28}H_{36}N_2O_3S$
- MW: 481
- [90134-00-2]

* 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. Nature, 355, 349(1992), [DOI: 10.1038/355349a0](#)
2. Biochim Biophys Acta, 1068(2), 149(1991), [DOI: 10.1016/0005-2736\(91\)90203-K](#)
3. PLoS ONE 8(10), e75678 (2013), [DOI: 10.1371/journal.pone.0075678](#)
4. Am J Phys Heart, 0(0), (2021), [DOI: 10.1152/ajpheart.00185.2021](#)

This datasheet was generated on January 21, 2026 at 05:03:30 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/di-4-anepps/>