Glowing products for science

DiA

DiA 4-(4-Dihexadecylaminostyryl)-N-methylpyridinium iodide) is a green fluorescent membrane dye which diffuses much faster than DiO in cell membranes. DiA and DiI (60010) have been used together for two color membrane staining.



Product Description

DiA 4-(4-Dihexadecylaminostyryl)-N-methylpyridinium iodide) is a green fluorescent membrane dye which diffuses much faster than DiO in cell membranes. DiA and DiI (60010) have been used together for two color membrane staining.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 491/613 nm
- Red solid soluble in DMSO
- Store at 4°C and protect from light, especially in solution
- C₄₆H₇₉IN₂
- MW: 787
- [114041-00-8]

References

1. Trends Neurosci 13, 14 (1990). 2. J Neurosci 12, 3494 (1992). 3. J Neurosci 15, 990 (1995).

This datasheet was generated on November 18, 2025 at 11:31:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dia-4-4-dihexadecylaminostyryl-n-methylpyridinium-iodide/

Product attributes CAS number Probe cellular localization

For live or fixed cells

114041-00-8

Membrane/cell surface, Membrane/vesicular

For fixed cells, For live/intact cells

Assay type/options

Co-cultures, Extended staining (several days to weeks)

Fixation options

Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining

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

Colors Green
Excitation/Emission 491/613 nm