

DiA

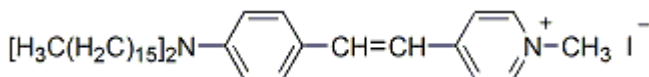
DiA 4-(4-Dihexadecylaminostyryl)-N-methylpyridinium iodide) is a green fluorescent membrane dye which diffuses much faster than DiO in cell membranes. DiA and Dil (60010) have been used together for two color membrane staining.



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 Dil ([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
- $C_{46}H_{79}IN_2$
- 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 January 12, 2026 at 07:40:42 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	114041-00-8
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	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
Colors	Green
Excitation/Emission	491/613 nm