



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 4-Di-1-ASP

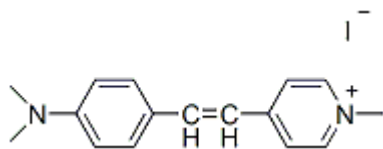
4-Di-1-ASP is a styryl dye that stains mitochondria of live cells. It has also been used to stain glioma cells in living brain tissues.



## Product Description

4-Di-1-ASP is a styryl dye that stains mitochondria of live cells. It has also been used to stain glioma cells in living brain tissues.

- $\lambda_{Ex}/\lambda_{Em} = 474/606$  nm (MeOH)
- Store at 4°C and protect from light, especially in solution
- $C_{16}H_{19}IN_2$
- MW: 366.2
- [68971-03-9]



## References

1. BioTechniques (2013) doi:10.2144/000113940
2. Journal of Biomolecular Screening (2016), doi:10.1177/1087057116646553

This datasheet was generated on June 9, 2026 at 07:47:12 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/4-4-dimethylamino-styryl-n-methylpyridinium-iodide-4-di-1-asp/>

## Product attributes

CAS number	68971-03-9
Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining
Cell permeability	Membrane permeant
Apoptosis/viability marker	Mitochondrial potential
Colors	Red
Excitation/Emission	474/606 nm