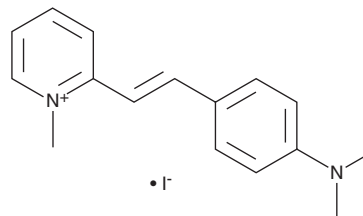


PRODUCT INFORMATION



2-DI-1-ASP Item No. 42328

CAS Registry No.: 2156-29-8
Formal Name: 2-[2-[4-(dimethylamino)phenyl]ethenyl]-1-methyl-pyridinium, monoiodide
Synonyms: DASPI, DASPMI
MF: C₁₆H₁₉N₂ • I
FW: 366.2
Purity: ≥98%
Ex./Em. Max: 474/605 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-DI-1-ASP is supplied as a solid. A stock solution may be made by dissolving the 2-DI-1-ASP in the solvent of choice, which should be purged with an inert gas. 2-DI-1-ASP is sparingly soluble (1-10 mg/ml) in DMSO.

Description

2-DI-1-ASP is a fluorescent mitochondrial probe.¹ It has been used to image actively respiring mitochondria *in vitro*. 2-DI-1-ASP displays excitation/emission maxima of 474/605 nm, respectively.²

References

1. Villa, A.M., and Doglia, S.M. Mitochondria in tumor cells studied by laser scanning confocal microscopy. *J. Biomed. Opt.* **9**(2), 385-394 (2004).
2. Swayne, T.C., Gay, A.C., and Pon, L.A. Visualization of mitochondria in budding yeast. *Methods Cell Biol.* **80**, 591-626 (2007).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/11/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM