

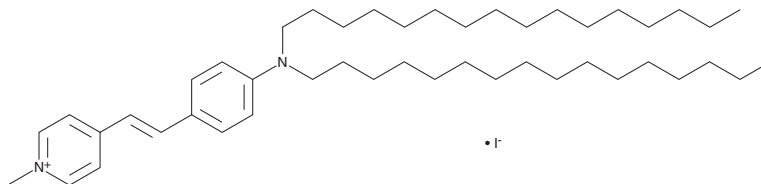
PRODUCT INFORMATION



DiA

Item No. 45034

CAS Registry No.: 114041-00-8
Formal Name: 4-[2-[4-(dihexadecylamino)phenyl]ethenyl]-1-methyl-pyridinium, monoiodide
Synonym: Di-16-ASP
MF: C₄₆H₇₉N₂ • I
FW: 787.1
Purity: ≥90%
Ex./Em.: 491/614 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

DiA is supplied as a solid. A stock solution may be made by dissolving the DiA in the solvent of choice, which should be purged with an inert gas. DiA is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1) and DMSO.

Description

DiA is a lipophilic, fluorescent dye.^{1,2} It displays excitation/emission maxima of 491/613 nm, respectively. DiA has been used in neuronal membrane tracing.¹

References

1. Ince-Dunn, G., Hall, B.J., Hu, S.C., *et al.* Regulation of thalamocortical patterning and synaptic maturation by NeuroD2. *Neuron* **49**(5), 683-695 (2006).
2. Neuhuber, W.L., Kressel, M., Stark, A., *et al.* Vagal efferent and afferent innervation of the rat esophagus as demonstrated by anterograde Dil and DiA tracing: Focus on myenteric ganglia. *J. Auton. Nerv. Syst.* **70**(1-2), 92-102 (1998).

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, 02/26/2026

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