

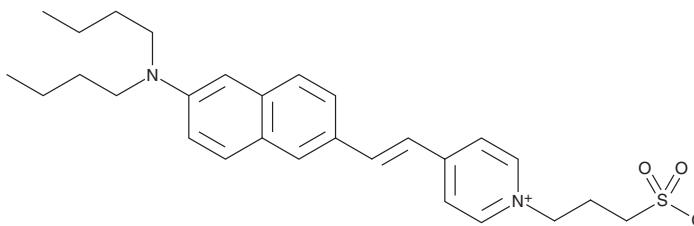
PRODUCT INFORMATION



Di-4-ANEPPS

Item No. 26801

CAS Registry No.: 90134-00-2
Formal Name: 4-[2-[6-(dibutylamino)-2-naphthalenyl]ethenyl]-1-(3-sulfopropyl)-pyridinium, inner salt
MF: C₂₈H₃₆N₂O₃S
FW: 480.7
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Di-4-ANEPPS is supplied as a solid. A stock solution may be made by dissolving the Di-4-ANEPPS in the solvent of choice. Di-4-ANEPPS is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

Di-4-ANEPPS is a potentiometric fluorescent dye that can be used to measure membrane potential.¹ It displays absorption/emission maxima of 495/705 nm, respectively.² The relative fluorescence changes proportionally to membrane potential at a rate of approximately 10% per 100 mV.¹ Di-4-ANEPPS has been used to measure membrane potential in a variety of tissue and cell types as well as artificial membranes.^{1,2}

References

1. Loew, L.M., Cohen, L.B., Dix, J., *et al.* A naphthyl analog of the aminostyryl pyridinium class of potentiometric membrane dyes shows consistent sensitivity in a variety of tissue, cell, and model membrane preparations. *J. Membr. Biol.* **130(1)**, 1-10 (1992).
2. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. John Wiley & Sons, Inc., Madison, NJ (2010).

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, 01/22/2019

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