

# Di-8-ANEPPS

## Membrane Potential Detector

Di-8-ANEPPS is a member of the ANEP class of membrane potential dyes. These dyes are weakly fluorescent in aqueous media, and become strongly fluorescent upon binding to lipophilic environments such as membranes. A change in the surrounding electronic field demonstrates a membrane potential-dependent shift in excitation spectra. Di-8-ANEPPS is less susceptible for cellular internalization than other ANEP dyes probably due to a sulfonate group. The response is sufficiently fast to detect transient (millisecond) potential changes in excitable cells, including single neurons, cardiac cells and intact brains. Wavelength Maxima: Excitation 498nm, Emission 713nm

Citations: 3

[View Online »](#)

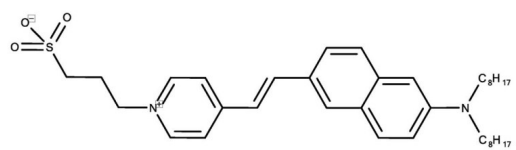
## Ordering Information

[Order Online »](#)

ENZ-52204	5mg
-----------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	4-[2-[6-(Diocetyl amino)-2-naphthalenyl]ethenyl]-1-(3-sulfopropyl)-pyridinium, inner salt
CAS	157134-53-7
Formula	$C_{36}H_{52}N_2O_3S$
MW	592.9
Purity	≥90% (HPLC)
Solubility	Soluble in DMSO.

Technical Info / Product Notes	This product is a member of the CELLESTIAL <sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL <sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.
--------------------------------	-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)