Dipicrylamine

Dipicrylamine (DPA) is a nonfluorescent anionic membrane dye. With its absorbance in the blue region of the visible spectrum, DPA can be used as a FRET quencher for a variety of fluorescent donor dyes.



Product attributes

CAS number	131-73-7
For live or fixed cells	For live/intact cells
Potential dependence	Slow response (translational) membrane potential dve

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Dipicrylamine (DPA) is a nonfluorescent anionic membrane dye. With its absorbance in the blue region of the visible spectrum, DPA can be used as a FRET quencher for a variety of fluorescent donor dyes. When applied to cells, DPA localizes at the interface of cytoplasmic membrane and aqueous phase in a manner sensitive to membrane potential change. A recent application of DPA is to measure cell membrane potential change by combining with the fluorescent donor dye DiOC16 (cat# 60038). The DiO/DPA system has been reported to produce a fluorescence signal change of >56% in HEK-293 cells and >25% in neuronal cultures and brain slices per 100 mV membrane potential change. Please also see the DiO/DPA Membrane Potential Detection Kit (30037)).

Yellow solid soluble in DMSO at >30mg/mL

- λabs (MeOH) = 406 nm
- Store at 4°C
- MWt: 439.21
- [131-73-7]

$$O_2N$$
 NO_2
 O_2N
 NO_2
 O_2N
 NO_2
 O_2N

References

1. J Neurosci 29(29), 9197 (2009).

This datasheet was generated on November 8, 2025 at 10:15:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dipicrylamine/