

## 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

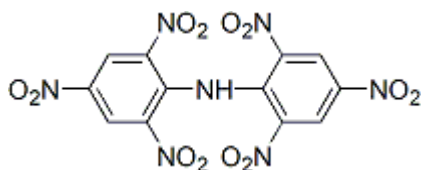
CAS number	131-73-7
For live or fixed cells	For live/intact cells
Potential dependence	Slow response (translational) membrane potential dye

## 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

- $\lambda_{\text{abs}}$  (MeOH) = 406 nm
- Store at 4 °C
- MWt: 439.21
- [131-73-7]



## References

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

This datasheet was generated on January 2, 2026 at 08:09:15 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dipicrylamine/>