

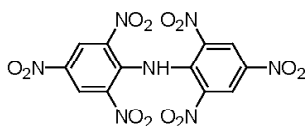
Product Information

Dipicrylamine

Catalog Number: 60037

Packaging Size: 25 mg

Molecular Information:

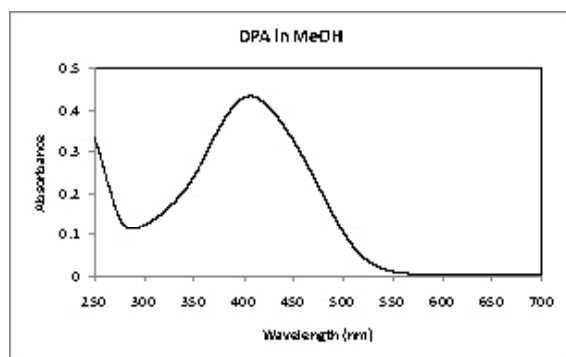


Mwt: 439.21; CAS#: [131-73-7]

Color and Form: Yellow solid

Solubility: Soluble in DMSO at >30mg/mL

Spectral Properties: λ_{abs} (MeOH) = 406 nm



Storage and Handling

Store at room temperature or below. Shelf-life is at least 1 year from the date of receipt.

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 (Bradley, et al. J. Neurosci. 29(29), 9197(2009)).

Reference:

1. Bradley, et al. J. Neurosci. 29(29), 9197(2009).
2. Chanda, et al. Nature 436, 852(2005).

Safety

See MSDS.

First aid

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Other related products

- 1) DiO/DPA membrane potential detection kit (cat# 30037).
- 2) DiOC16(3) Cat# 60038).
- 3) Di-4-ANEPPS (Cat# 61010)
- 4) SynpatoGreen C4 (or FM1-43) (Cat# 70020)
- 5) Tetrodotoxin (Cat# 00061, 00062)
- 6) Fluo-3, AM (Cat# 50014)

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.