

Product datasheet

DAPI Staining Solution ab228549

★★★★★ [1 Abreviews](#) [166 References](#) [2 Images](#)

Overview

Product name DAPI Staining Solution

Sample type Adherent cells, Suspension cells

Product overview DAPI Staining Solution (ab228549) is a fluorescent stain for labeling DNA in fluorescence microscopy. Since DAPI passes through an intact cell membrane, it can be used to stain live cells and fixed cells.

Molecular Weight: 350.25

Notes DAPI's absorption maximum when bound to double-stranded DNA is at 358 nm and its emission maximum is at 461 nm. Although is not as strongly fluorescent as when bound to DNA, it also binds to RNA, shifting its emission to around 500 nm.

DAPI's blue emission is convenient for multiplexing assays since there is very little fluorescence overlap between DAPI and green-fluorescent molecules like FITC and green fluorescent protein (GFP), or red-fluorescent stains like Texas Red.

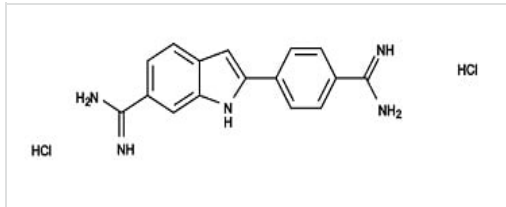
Platform Fluorescence microscope

Properties

Storage instructions Store at -20°C. Please refer to protocols.

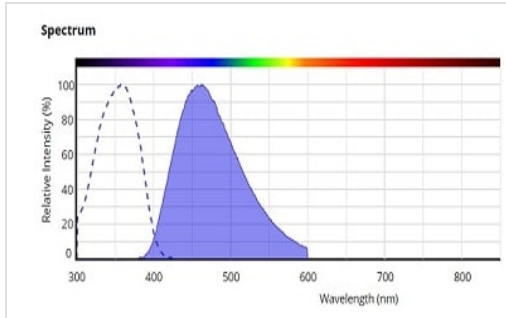
Components	2 ml
DAPI (10 mM solution in water)	1 x 2ml

Images



Chemical structure of DAPI.

Chemical structure of DAPI.



Spectrum of DAPI.

Spectrum of DAPI.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors