

## ab228549 - DAPI Staining Solution

For labeling DNA in fluorescence microscopy.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/ab228549](http://www.abcam.com/ab228549)  
(use [www.abcam.cn/ab228549](http://www.abcam.cn/ab228549) for China, or [www.abcam.co.jp/ab228549](http://www.abcam.co.jp/ab228549) for Japan)

### General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

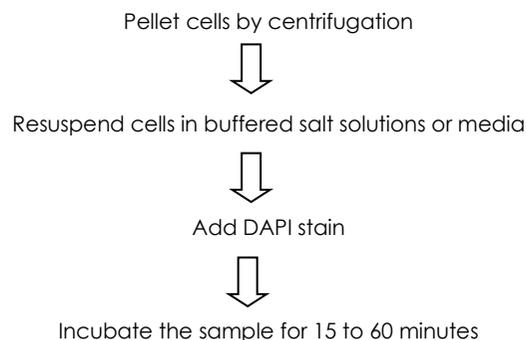
For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:  
[www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines)

For typical data produced using the assay, please see the assay kit datasheet on our website.

### 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.

### Assay Summary



### Materials Supplied

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
DAPI [4,6-Diamidino-2-phenylindole, dihydrochloride] (10 mM solution in water)	2 mL	-20°C	-20°C

### Storage and Stability

Store kit at -20°C and desiccated in the dark immediately on receipt and check below for storage for individual components. Kit can be stored for 6 months from receipt, if components have not been reconstituted.

### Materials Required, Not Supplied

These materials are not included, but will be required to successfully perform this assay:

- Buffered salt solutions or media, with optimal dye binding at pH 7.4.

### Reagent Preparation

Briefly centrifuge small vial at low speed prior to opening.

### **DAPI 4,6-Diamidino-2-phenylindole, dihydrochloride] (10 mM solution in water)**

Ready to use as supplied.

### Assay procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
- The following procedure can be adapted for most cell types. Growth medium, cell density, the presence of other cell types and other factors may influence staining.
- Residual detergent on glassware may also affect real or apparent staining of many organisms, causing brightly stained material to appear in solutions with or without cells present.

DAPI staining is normally performed after all other staining.

1. Pellet cells by centrifugation and resuspend the cells in buffered salt solutions or media, with optimal dye binding at pH 7.4.
2. Adherent cells in culture may be stained *in situ* on cover slips or in the cell culture wells.
3. Add DAPI stain using the concentrations between 0.5 and 5  $\mu$ M and incubate it for 15 to 60 minutes as a guide.
4. In initial experiments, it may be best to try several dye concentrations over the entire suggested range to determine the concentration that yields optimal staining.

### FAQs / Troubleshooting

General troubleshooting points can be found at [www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines).

### Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Alexa Fluor® is a registered trademark of Life Technologies.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)