

ab228551 Hoechst 33342 Staining Dye Solution

For labeling DNA in fluorescence microscopy. This product is for research use only and is not intended for diagnostic use.

Overview

Hoechst 33342 Staining Dye Solution (ab228551) is a fluorescent stain for labeling DNA in fluorescence microscopy. This product may be used in fluorescence microscopy, microplate, cuvette, and flow cytometry applications. It can also be used to detect the contents of a sample DNA by plotting a standard emission-to content curve.

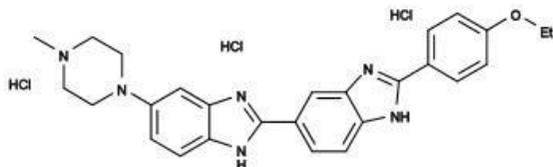


Figure 1. Chemical structure of Hoechst 33342.

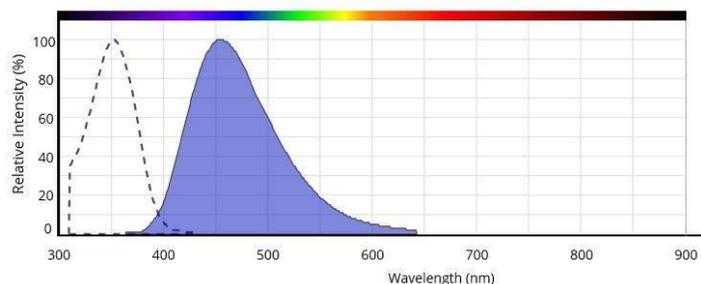


Figure 2. Spectrum of Hoechst 33342.

Materials Supplied and Storage

Store kit at -20°C in the dark immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Item	Quantity	Storage temperature (before prep)
Hoechst 33342 (20 mM Solution in Water)	5 mL	-20°C

Materials Required, Not Supplied

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

1. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

Hoechst 33342 (20 mM Solution in Water): Ready to use as supplied.

2. 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.

2.1 Pellet cells by centrifugation.

2.2 Resuspend the cells in buffered salt solutions or media, with optimal dye binding at pH 7.4.

2.3 Adherent cells in culture may be stained *in situ* on cover slips or in the cell culture wells.

2.4 Add Hoechst stain using the concentrations between 0.5 and 5 μ M and incubate it for 15 to 60 minutes as a guide.

2.5 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.

General troubleshooting points can be found at www.abcam.com/assaykitguidelines.

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)