

Product datasheet

Hoechst 33342 Staining Dye Solution ab228551

★★★★★ [1 Abreviews](#) [16 References](#) [2 Images](#)

Overview

Product name Hoechst 33342 Staining Dye Solution

Sample type Adherent cells, Suspension cells

Product 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-tocontent curve.

Molecular Weight: 561.93

Spectral Properties: Excitation = 350 nm; Fluorescence = 461 nm.

Notes The Hoechst stains are a family of fluorescent stains for labeling DNA in fluorescence microscopy. Because these fluorescent stains label DNA, they are also commonly used to visualize nuclei and mitochondria. Two of these closely related bis-benzimides are commonly used: Hoechst 33258 and Hoechst 33342. Both dyes are excited by ultraviolet light at around 350 nm, and both emit blue/cyan fluorescence light with an emission maximum at 461 nm. The Hoechst stains may be used on live or fixed cells, and are often used as a substitute for another nucleic acid stain, DAPI. The key difference between them is that the additional ethyl group of Hoechst 33342 makes it more lipophilic, and thus more readily to cross intact cell membranes. In some applications, Hoechst 33258 is significantly less permeant than Hoechst 33342.

Tested applications **Suitable for:** FM, Flow Cyt

Properties

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

Components	5 ml
Hoechst 33342 (20 mM Solution in Water)	1 x 5ml

Applications

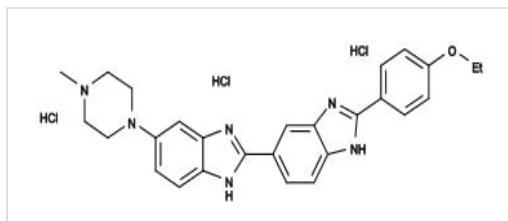
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab228551 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

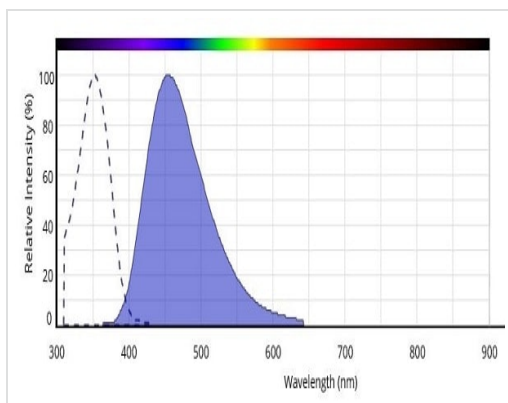
Application	Abreviews	Notes
FM		Use at an assay dependent concentration.
Flow Cyt		Use at an assay dependent concentration.

Images



Chemical structure of Hoechst 33342.

Chemical structure of Hoechst 33342.



Spectrum of Hoechst 33342.

Spectrum of Hoechst 33342.

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