Hoechst Stains

Table 1. Contents and storage information.

Product	Amount	Concentration	Storage	Stability
Hoechst 33258, pentahydrate (bis-benzimide) FluoroPure™ grade (MW 623.96) (H21491)	100 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.
Hoechst 33258, pentahydrate (bis-benzimide) (MW 623.96) (H1398)	100 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.
Hoechst 33258, pentahydrate (bis-benzimide) (MW 623.96) (H3569)	10 mL	10 mg/mL (16.0 mM) solution in water	Store at 2–6°C Protect from light	When stored as directed, product is stable for 6 months.
Hoechst 33342, trihydrochloride, trihydrate, FluoroPure™ grade (MW 615.99) (H21492)	100 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.
Hoechst 33342, trihydrochloride, trihydrate, (MW 615.99) (H1399)	100 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.
Hoechst 33342, trihydrochloride, trihydrate, (MW 615.99) (H3570)	10 mL	10 mg/mL (16.2 mM) solution in water	Store at 2–6°C Protect from light	When stored as directed, product is stable for 6 months.
Hoechst 34580, (MW 560.96) (H21486)	5 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.
nuclear yellow (Hoechst S769121, trihydrochloride, trihydrate, (MW 651.01) (N21485)	10 mg	NA	• ≤25°C • Protect from light	When stored as directed, product is stable for at least 12 months.

Approximate Fluorescence Excitation and Emission, in nm: Excitation/Emission bound to DNA: Hoechst 33258 352/461 nm (see Figure 2); Hoechst 33342 350/461 nm; Hoechst 34580 392/440 nm (see Figure 3); Nuclear Yellow 355/495 nm.

Introduction

The blue fluorescent Hoechst dyes (Figure 1) are cell permeable nucleic acid stains that have multiple applications, including sensitive detection (>3 ng) of DNA in the presence of RNA in agarose gels, 1 automated DNA determination, 2 sensitive determination of cell number 3,4 and chromosome sorting.⁵ The fluorescence of these dyes is very sensitive to DNA conformation and chromatin state in cells. Consequently, they can detect gradations of nuclear damage. The Hoechst dyes are useful vital stains for the flow cytometric recognition of DNA damage ^{6,7} and other viability measurements by monitoring the emission spectral shifts of the dyes.⁸ These bisbenzimidazole derivatives are supravital minor groove-binding DNA stains with AT selectivity. The dyes

bind to all nucleic acids, but AT-rich dsDNA strands enhance fluorescence ~2-fold greater than GC-rich strands. This property has been used to identify Q-bands in chromosomes (Q-bands: AT-rich chromosome regions that fluoresce brightly when stained with the dye quinacrine). 10

$$\begin{array}{c} \text{CH}_{3}\text{NH} \\ \text{N} \\ \text{N} \\ \text{H} \\$$

Figure 1. Structure of Hoechst dyes.

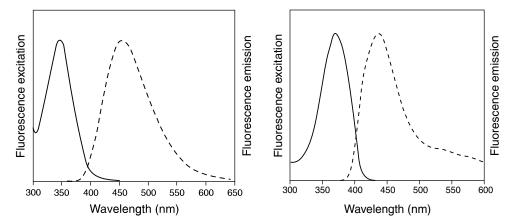


Figure 2. Fluorescence excitation and emission spectra of Hoechst 33258 bound to DNA

Figure 3. Fluorescence excitation and emission spectra of Hoechst 34580 bound to DNA.

BrdU Quenching

Hoechst 33342 and 33258 are quenched by BrdU. 11,12 Fluorescence is enhanced upon binding to dsDNA at stretches of at least three AT base pairs, but no binding to stretches of two or more GC base pairs. The Hoechst dyes requires a [dA-dT]₃-[dG-dC]₁ sequence to enhance fluorescence, with binding to the bottom of the minor groove as a prerequisite. It is postulated that the bromine on BrdU (or the chlorine on CldU) deforms the minor groove, such that the dye molecule can no longer reach its optimal binding site. It has shown that the Hoechst dyes still bind (with an even higher affinity) to BrdU substituted DNA, but no fluorescence ensues. 13 This has been used extensively to monitor cell cycle progression.

Hoechst 33258 is slightly more water soluble than Hoechst 33342, but both have been used extensively to stain live cells. The products may be used in fluorescence microscopy, microplate, cuvette, and flow cytometry applications. Nuclear yellow (Hoechst S769121) is more commonly used as a neuronal retrograde tracer. 14-16

Preparing Stock Solutions of Hoechst Dyes

The solid dyes may be dissolved in either water, dimethylformamide (DMF), or DMSO to make concentrated stock solutions up to 10 mg/mL. Stock solutions may be stored refrigerated or frozen, protected from light. Note: The Hoechst stains should not be resolubilized in phosphate-buffered saline (PBS), but dilute solutions of the dye may be used with PBS or other phosphate-containing buffers. Solutions of Hoeschst dye should be stored at 2–6°C, protected from light. Stock solutions in water are stable for at least 6 months when refrigerated. For long-term storage the stock solution can be aliquoted and stored at $\leq -20^{\circ}$ C.

Caution

The Hoechst strains are known mutagens and should be handled with care. The dye must be disposed of safely and in accordance with applicable regulations.

Fluorescence Spectral **Characteristics**

The blue fluorescent Hoechst 33258, 33342, 34580 and nuclear yellow dyes may be efficiently excited with a xenon or mercury-arc lamp or with a UV laser; Hoechst 34580 may also be efficiently excited with non-UV wavelengths (>360 nm) such as the 405 nm laser line. They may all be detected using the common DAPI filter, blue GFP filters or the Semrock BrightLine® Alexa Fluor® 350 filter set.

Basic Protocol for Staining Cells

The following procedure can be adapted for most cell types. Note that different concentration ranges for the Hoechst dyes are suggested depending on the cell type (see Table 2). 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. Glassware should be washed in a mild detergent and rinsed with hot tap water followed by several rinses with deionized, distilled water.

Pellet cells by centrifugation and resuspend in buffered salt solutions or media, with optimal dye binding at pH 7.4. Adherent cells in culture may be stained *in situ* on coverslips. Add Hoechst stain using the concentrations listed in Table 2 as a guide. 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. Unbound dye has its maximum fluorescence emission in the 510-540 nm range, this green fluorescence may be observed on samples using too high a concentration of dye.

Table 2. Recommended	conditions for staining	g cells with Hoechst stains.

Cell Type	Hoechst Dye Concentration	Incubation Conditions
Bacteria	0.1 to 12 μg/mL	10 to 30 minutes
Live animal cells	0.2 to 5 μg/mL	20 to 30 minutes
Fixed animal cells	0.2 to 2 μg/mL	1 to 15 minutes

References

^{1.} Nucleic Acids Res 15, 10589 (1987); 2. Anal Biochem 147, 462 (1985); 3. Anal Biochem 131, 538 (1983); 4. J Histochem Cytochem 29, 326 (1981); 5. Cytometry 3, 145 (1982); 6. Cytometry 11, 386 (1990); 7. Cancer Res. 48, 5742 (1988); 8. Cytometry 11, 239 (1990); 9. BBA 949, 158 (1988); 10. Chromosoma 46, 255 (1974); 11. J Histochem Cytochem 24, 24 (1976); 12. J Histochem Cytochem 25, 913 (1977); 13. J Mol Biol 315, 1049 (2002); 14. Neurosci Lett 18, 19 (1980); 15. Exp Brain Res 40, 383 (1980); 16. J Histochem Cytochem 30, 123 (1982).

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat #	Product Name	Unit Size
F2962	FluoReporter® Blue Fluorometric dsDNA Quantitation Kit *200-2000 assays*	1 kit
N21485	nuclear yellow (Hoechst S769121, trihydrochloride, trihydrate)	10 mg
H1398	Hoechst 33258, pentahydrate (bis-benzimide)	100 mg
H3569	Hoechst 33258, pentahydrate (bis-benzimide) *10 mg/mL solution in water*	10 mL
H21491	Hoechst 33258, pentahydrate (bis-benzimide) *FluoroPure™ grade*	100 mg
H1399	Hoechst 33342, trihydrochloride, trihydrate	100 mg
H3570	Hoechst 33342, trihydrochloride, trihydrate *10 mg/mL solution in water*	10 mL
H21492	Hoechst 33342, trihydrochloride, trihydrate *FluoroPure™ grade*	100 mg
H21486	Hoechst 34580	5 mg
V13244	Vybrant® Apoptosis Assay Kit #5 *Hoechst 33342/propidium iodide* *200 assays*	1 kit
V23201	Vybrant® Apoptosis Assay Kit #7 *Hoechst 33342/YO-PRO®-1/propidium iodide* *200 assays*	1 kit

Contact Information

Molecular Probes, Inc.

29851 Willow Creek Road Eugene, OR 97402 Phone: (541) 465-8300 Fax: (541) 335-0504

Customer Service:

6:00 am to 4:30 pm (Pacific Time) Phone: (541) 335-0338 Fax: (541) 335-0305 probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209 Order Fax: (800) 438-0228

Technical Service:

8:00 am to 4:00 pm (Pacific Time) Phone (541) 335-0353 Toll-Free (800) 438-2209 Fax: (541) 335-0238 probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd. 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Phone: +44 (0) 141 814 6100 Fax: +44 (0) 141 814 6260 Email: euroinfo@invitrogen.com Technical Services: eurotech@invitrogen.com

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Assistance Department in Eugene, Oregon.

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License

For research use only. Not intended for any animal or human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402. Tel: (541) 465-8300. Fax: (541) 335-0504.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ° are registered with the U.S. Patent and Trademark Office.

Copyright 2005, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.