

Revised: February 21, 2019

# **Product Information**

# Hoechst 33258

| Catalog number | Product                           | MW     | Size   |
|----------------|-----------------------------------|--------|--------|
| 40044          | Hoechst 33258, 10 mg/mL in $H_2O$ | 533.88 | 10 mL  |
| 40045          | Hoechst 33258, pentahydrate       | 623.96 | 100 mg |

# Storage and Handling

Store Hoechst 33258, 10 mg/mL in  $H_2O$  at 4°C, protected from light. Store Hoechst 33258, pentahydrate solid desiccated at -20°C, protected from light. Product is stable for at least one year from date of receipt when stored as recommended. It is not recommended to store working solutions of Hoechst dye, because the dye will be lost to precipitation or adsorption to the container over time.

# **Molecular Information**

C25H27CI3N6O

## CAS number: 23491-45-4

Color and form:

40044: Pale yellow solution 40045: Yellow solid

40045: Yellow solid

Solubility: Soluble in water up to 20 mg/mL

Absorption/Emission: 352/461 nm (with DNA) (Figure 2)



Figure 1. Hoechst 33258, pentahydrate.

# **Product Description**

Hoechst dyes are widely-used cell membrane-permeant blue fluorescent DNA binding dyes that can be used for nuclear staining of live or fixed cells. The dyes have minimal fluorescence in solution, but become brightly fluorescent upon binding to DNA. Therefore, they can be used to stain cells without a wash step. The staining is very stable and non-toxic to live cells for several days or longer.

Hoechst 33342 and Hoechst 33258 are structurally similar minor-groove DNA binding dyes that perform comparably as nuclear counterstains. Hoechst 33528 is slightly more water soluble than Hoechst 33342, but both dyes are highly cell membrane permeant and widely used in cell cycle studies and as nuclear counterstains for live or fixed cells.

## References

1) Science 282, 296 (1998); 2) Biotech. Histochem. 69, 152 (1994); 3) Exp. Cell Res. 211, 322 (1994).





Figure 2. Normalized excitation and emission spectra of Hoechst 33258 with DNA.

# **Staining Protocols**

## Live cell staining

Below we provide two protocols for staining live cells with Hoechst. Staining by medium exchange results in uniform exposure of cells to probe. However, for some cell types, morphology or viability may be affected by medium exchange. In addition, floating dead cells may be lost during medium removal, and suspension cells must be collected by centrifugation to exchange the medium. Direct addition of 10X probe is a convenient staining method that doesn't require medium exchange, but care must be taken to mix immediately yet gently to avoid high transient probe concentration or disruption of cells by pipetting. Note that we do not recommend adding highly concentrated dye directly to cells in culture, as this will result in local areas of high dye exposure.

## Live cell staining by medium exchange

- 1. Dilute Hoechst dye to 1 ug/mL in fresh, complete cell culture medium. Hoechst can be combined with other fluorescent probes.
- 2. Remove medium from the cells and replace with fresh medium containing Hoechst dye.
- Incubate cells at room temperature or 37°C for 5-15 minutes, then image. Note: Washing is not necessary for specific staining, but nuclear staining is stable after washing.

## Live cell staining by direct addition of 10X probe

- Prepare 10X dye solution by diluting Hoechst dye to 10 ug/mL in fresh, complete culture medium. Hoechst can be combined with other fluorescent probes, which should be diluted to 10 times the final desired concentration.
- 2. Without removing the medium from the cells, add 1/10 volume of 10X dye solution directly to the well.
- Immediate mix thoroughly by gently pipetting the medium up and down. For larger well sizes (e.g., 24-well to 6-well plates), the plate can be gently swirled to mix.
- Incubate cells at room temperature or 37°C for 5-15 minutes, then image. Note: Washing is not necessary for specific staining, but nuclear staining is stable after washing.

## Staining of fixed cells or tissue sections

- Add Hoechst dye to PBS at 1 ug/mL. Hoechst can be included together with antibodies or other probes, and can be diluted in buffers with detergent or blocking agents if convenient.
- Add the PBS with dye to cells or tissue sections and incubate at room temperature for at least 5 minutes.
- Image the samples; washing is optional but not required. Note: Samples can be stored at 4°C after staining and before imaging.

## Staining bacteria or yeast

Hoechst dyes stain bacteria more dimly than mammalian cells. Live or killed bacteria can be stained with 12-15 ug/mL Hoechst dye in PBS or 150 mM NaCl for 30 minutes at room temperature. The dyes tend to stain dead cells more brightly than live cells.

In *S. cerevisiae*, Hoechst dyes stain both the nucleus and cytoplasm. When used to stain yeast at 12-15 ug/mL in PBS, Hoechst preferentially stains dead cells, with dim staining of live cells.

# **Related Products**

| Catalog<br>number | Product  |
|-------------------|--|
| 40046             | Hoechst 33342, 10 mg/mL in dH2O                          |
| 40047             | Hoechst 33342, trihydrochloride trihydrate               |
| 40083             | NucSpot® 470   |
| 40081             | NucSpot® Live 488  |
| 40082             | NucSpot® Live 650  |
| 40085             | NucSpot® Far-Red   |
| 40060             | RedDot™1 Far-Red Nuclear Stain                           |
| 40061             | RedDot™2 Far-Red Nuclear Stain                           |
| 40084             | 7-AAD Solution, 1 mg/mL                                  |
| 40048             | Propidium lodide Buffer, 50 ug/mL                        |
| 40009             | DAPI, dilactate  |
| 40011             | DAPI, dihydrochloride                                    |
| 40043             | DAPI, dilactate 10 mg/mL in H <sub>2</sub> O             |
| 23001             | EverBrite™ Mounting Medium                               |
| 23002             | EverBrite™ Mounting Medium with DAPI                     |
| 23003             | EverBrite™ Hardset Mounting Medium                       |
| 23004             | EverBrite <sup>™</sup> Hardset Mounting Medium with DAPI |
| 23008             | Drop-n-Stain EverBrite™ Mounting Medium                  |
| 23009             | Drop-n-Stain EverBrite™ Mounting Medium with DAPI        |
| 30068             | ViaFluor® 405 SE Cell Proliferation Kit                  |
| 30086             | ViaFluor® 488 SE Cell Proliferation Kit                  |
| 70065             | LipidSpot™ 488 Lipid Droplet Stain                       |
| 70069             | LipidSpot™ 610 Lipid Droplet Stain                       |
| 70066             | LysoView™ 405 Lysosome Stain                             |
| 70067             | LysoView™ 488 Lysosome Stain                             |
| 70061             | LysoView™ 540 Lysosome Stain                             |
| 70058             | LysoView™ 633 Lysosome Stain                             |
| 70059             | LysoView™ 650 Lysosome Stain                             |
| 70070             | MitoView™ 405 Mitochondrial Stain                        |
| 70054             | MitoView™ Green Mitochondrial Stain                      |
| 70055             | MitoView™ 633 Mitochondrial Stain                        |
| 70075             | MitoView™ 650 Mitochondrial Stain                        |
| 70068             | MitoView™ 720 Mitochondrial Stain                        |
| 70064             | ViaFluor® 405 Live Cell Microtubule Stain                |
| 70062             | ViaFluor® 488 Live Cell Microtubule Stain                |
| 70063             | ViaFluor® 647 Live Cell Microtubule Stain                |
| 30090             | CellBrite™ Fix 488 Membrane Stain                        |
| 30088             | CellBrite™ Fix 555 Membrane Stain                        |
| 30089             | CellBrite™ Fix 640 Membrane Stain                        |
| 30092-<br>30099   | MemBrite™ Fix Cell Surface Staining Kits                 |
| 22023             | Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative       |
| 22020             | 10X Phosphate-Buffered Saline (PBS)                      |

Please visit our website at www.biotium.com for information on our life science research products, including novel live cell stains for nuclei, mitochondria, and other organelles, fluorescent CF® dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.