

Hoechst 33342-Biotin Conjugate

Catalog Number: 17675

Unit Size: 1 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

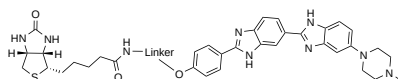
Chemical Properties

Appearance Solid

Molecular Weight 1307.61

Soluble In Water

Chemical Structure



Spectral Properties

Excitation Wavelength 352 nm

Emission Wavelength 454 nm

Applications

Hoechst 33342-Biotin Conjugate is a fluorescent nucleic acid stain that combines the DNA-binding properties of Hoechst 33342 with a biotin functional group for downstream detection. Hoechst dyes bind to the minor groove of double-stranded DNA, with preference for AT-rich regions, and are commonly used for nuclear staining in live or fixed cells due to their cell-permeable nature.

The biotin conjugation enables interaction with streptavidin or avidin-based detection systems, allowing signal amplification or probe localization in biochemical and imaging applications. This conjugate can be used in workflows that combine fluorescence imaging with affinity-based detection, including microscopy, flow cytometry, and nucleic acid-related assays.