

Helixyte™ Green Maleimide

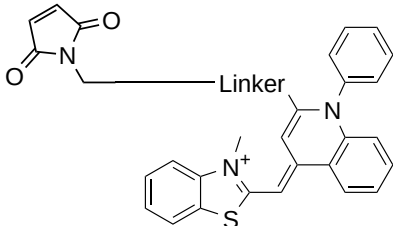
Catalog Number: 17614

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid yellow
Molecular Weight	973.05
Soluble In	Water
Chemical Structure	

Spectral Properties

Excitation Wavelength	498 nm
Emission Wavelength	522 nm

Applications

Helixyte™ Green is an excellent nucleic acid sensor that exhibits large fluorescence enhancement upon binding to dsDNA. It has the same spectral and DNA-binding properties as SYBR Green® (SYBR Green® is the trademark of ThermoFisher). AAT Bioquest's scientists developed Helixyte™ Green maleimide to enable the development of site-specific DNA probes. Helixyte™ Green maleimide is an excellent building block that readily reacts with thiol-containing biological molecules such as peptides, antibodies and thiol-modified oligos. Thus, it can be used to develop a variety of site-directed DNA probes. Helixyte™ Green is an excellent fluorogenic probe for quantifying DNAs with greatly improved sensitivity and selectivity. Helixyte™ Green is an ultra-sensitive fluorescent nucleic acid stain for quantitating double-stranded DNA (dsDNA) in molecular biological procedures such as cDNA synthesis for library production and DNA fragment purification for subcloning, as well as diagnostic applications, such as quantitating DNA amplification products and primer extension assays.