

# Helixyte™ Green NHS ester

Catalog Number: 17612

Unit Size: 1 mg

## Product Details

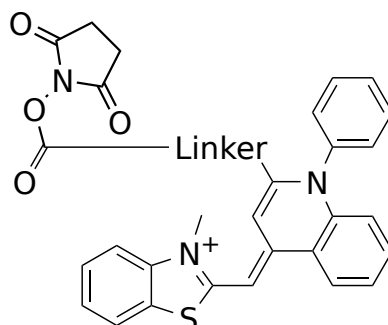
---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Solid red
Molecular Weight	947.99
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 NHS ester to enable the development of site-specific DNA probes. Helixyte™ Green NHS ester readily reacts with amino-containing biological molecules such as peptides, antibodies and amino-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. Using the Helixyte™ Green dsDNA Quantifying Reagent, you can selectively detect as little as 25 pg/ml of dsDNA in the presence of ssDNA, RNA, and free nucleotides.