

# Helixyte™ Green Amine

Catalog Number: 17613

Unit Size: 1 mg

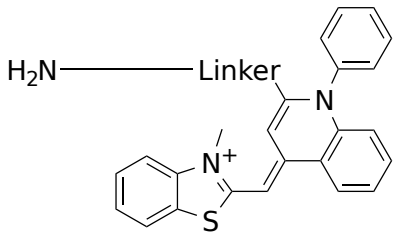
## Product Details

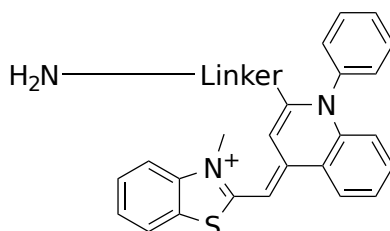
---

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

## Chemical Properties

---

Appearance	Solid
Molecular Weight	1007.03
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 amine to enable the development of site-specific DNA probes. Helixyte™ Green amine is an excellent building block that readily reacts with carbonyl-containing biological molecules such as peptides, antibodies and other proteins. 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.