

## Product Information Sheet

### Ordering Information

Product Number: 251  
Product Name: Helix Fluor™ 575, NHS ester  
Unit Size: 1 mg  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical, Physical and Spectral Properties

Molecular Weight: 770.31  
Appearance: Dark red solid  
Soluble in: DMSO  
Excitation Wavelength: 546 nm  
Emission Wavelength: 575 nm

### Application Notes

FAM, TET, VIC, JOE, NED, TAMRA and ROX fluorescent dyes are frequently used for labeling oligos. These fluorescent oligo conjugates are often used for nucleic acid sequencing and nucleic acid amplification. Our Helix Fluor™ 575 is an excellent alternative to the NED dye for preparing fluorescent oligos.