

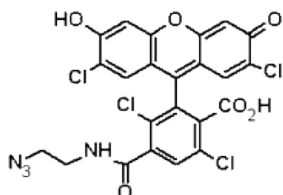
Product Information Sheet

Ordering Information

Product Number: 244
Product Name: 6-TET Azide
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 582.18
Appearance: Red solid
Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 521 nm
Emission Wavelength: 536 nm

Application Notes

6-TET is widely used in nucleic acid sequencing and related research. 6-TET azide provides a convenient method for labeling alkyne-modified oligos through the well-known click chemistry.