

Product Information Sheet

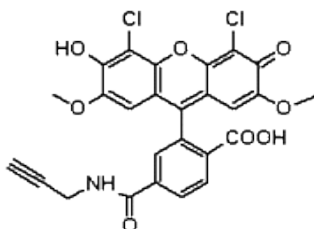
Ordering Information

Product Number: 249
Product Name: 6-JOE alkyne
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 542.32

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 520 nm
Emission Wavelength: 548 nm

Application Notes

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