

BG-TMR Conjugate

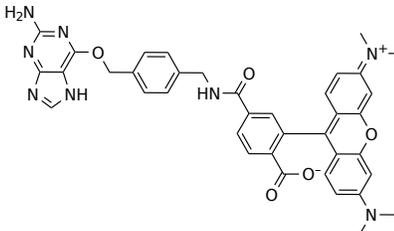
Catalog Number: 12558

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid red purple
Molecular Weight	682.74
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Site-specific labeling of cellular proteins with chemical probes is a powerful tool for studying protein functions in living cells. The SNAP-tag is a modified form of the DNA repair enzyme, human O6-alkylguanine-DNA-alkyltransferase (AGT) and undergoes a self-labeling reaction to form a covalent bond with O6-benzylguanine (BG) derivatives. BG-TMR conjugate is a fluorescent SNAP substrate that can be used to label SNAP-tag-fused target proteins. In the labeling reaction, the substituted benzyl group of the substrate is covalently attached to the SNAP-tag. SNAP-TMR substrate enables the specific, covalent attachment of TMR fluorophore to a protein of interest. The labeled proteins can be imaged with TRITC or Cy3 filter set.