

# Psoralen TMP Alkyne

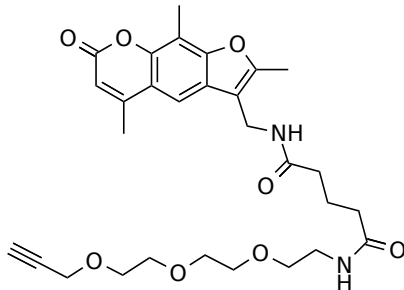
Catalog Number: 39063

Unit Size: 1 mg

## Product Details

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

## Chemical Properties

Appearance	Solid white
Molecular Weight	540.61
Soluble In	DMSO
Chemical Structure	

## Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

Psoralen TMP Alkyne is an excellent building block for developing site-specific nucleic acid probes via the well-known click chemistry, i.e., Copper-Catalyzed Azide-Alkyne Cycloaddition (CuAAC). It can be readily conjugated to azido-containing biomolecules (such as azido-modified antibodies, peptide azides and azido-modified oligos). Psoralens and their derivatives (such as 8-TMP and 4,5',8-TMP) are well known to have unique crosslinking features to DNA. Psoralen TMP Alkyne is an azide-reactive TMP derivative. It is an excellent building block for preparing TMP-labeled oligos from an azido-modified oligo. The TMP-conjugated oligonucleotides can be used for sequence-specific crosslinking with a target DNA, thus enabling the application of psoralen-conjugated molecules in gene transcription inhibition, gene knockout, and other genomic applications.