

## Psoralen MOP Alkyne

Catalog Number: 39058

Unit Size: 1 mg

### Product Details

---

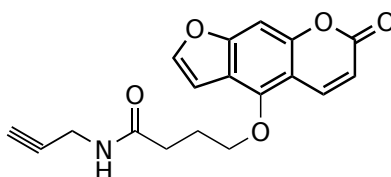
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	325.32
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

Psoralen MOP 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-MOP and 4,5'8-TMP) are well known to have unique crosslinking features to DNA. However, psoralen monomers do not have sequence-specific crosslinking ability with a target DNA. Nakao et al found that 5-MOP more effectively crosslinks DNA compared to the well-known 8-MOP (<https://doi.org/10.1002/cbic.202200789>). Psoralen MOP Alkyne is an azide-reactive 5-MOP derivative. It is an excellent building block for preparing MOP-labeled oligos from an azido-modified oligo. The 5-MOP-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.