

Psoralen TMP Succinimidyl Ester

Catalog Number: 39060

Unit Size: 5 mg

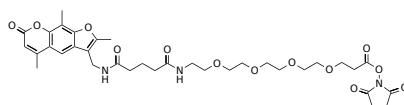
Product Details

| | |
|--------------------|--|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date | 12 months upon receiving |

Chemical Properties

| | |
|------------------|-------------|
| Appearance | Solid white |
| Molecular Weight | 715.74 |
| Soluble In | DMSO |

Chemical Structure



Spectral Properties

| | |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength | N/A |

Applications

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 Succinimidyl Ester is an amino-reactive TMP derivative. It is an excellent building block for preparing TMP-labeled oligos from the readily available amino-modified oligos. 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. Psoralen TMP Succinimidyl Ester may also be used for preparing site-specific DNA/RNA probes via the conjugations with amino-containing biomolecules such as antibodies.