

BG Succinimidyl Ester

Catalog Number: 12564

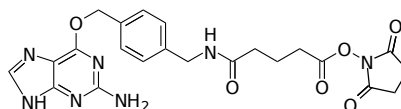
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid light yellow
Molecular Weight	481.47
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

SNAP protein labeling system enables the specific, covalent attachment of virtually any molecule to a protein of interest. The SNAP-tag is a protein based on human O6-alkylguanine-DNA-alkyltransferase (hAGT). SNAP-tag substrates are fluorophores, biotin or beads conjugated to guanine or chloropyrimidine leaving groups via a benzyl linker. In the labeling reaction, the substituted benzyl group of the substrate is covalently attached to the SNAP-tag. BG Succinimidyl Ester is an amine-reactive building block for preparing of SNAP substrates from amine-containing labels or biological targets. BG Succinimidyl Ester building block facilitates the development of various SNAP-tag substrates for labeling SNAP-tag fusion proteins for a wide range of biological applications.