

BG-Malaimide

Catalog number: 12566

Unit size: 1 mg

Product Details

Storage Conditions Freeze (<math>-15\text{ }^{\circ}\text{C}</math>), Minimize light exposure

Expiration Date 24 months upon receiving

Chemical Properties

Appearance Solid

Soluble In DMSO

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 Maleimide is a thiol-reactive building block for preparing of SNAP substrates from thiol-containing labels or biological targets. BG Maleimide building block facilitates the development of various SNAP-tag substrates for labeling SNAP-tag fusion proteins for a wide range of biological applications.