

DSG [Di(N-succinimidyl) glutarate] *CAS 79642-50-5*

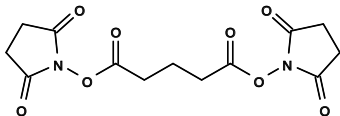
Catalog Number: 4520

Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid off-white
Molecular Weight	326.26
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

DSG is disuccinimidyl glutarate, a simple homobifunctional crosslinker based on the amine-reactive N-hydroxysuccinimide (NHS) ester group at both ends of a 5-atom (7.7 angstrom) spacer arm. It is used for crosslinking the molecules that have amine moiety. DSG is cell membrane-permeable. It is used in ChIP assays to determine DNA-protein binding interactions. DSG crosslinker must be dissolved in an organic solvent, such as DMSO or DMF, then added to an aqueous crosslinking reaction. Our in-house research demonstrated the DSG-crosslinked peptide conjugates have much better stability and tend to generate the antibodies of high specificity.