

EDTA Succinimidyl Ester

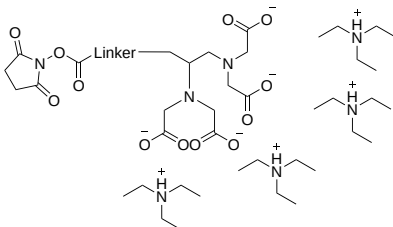
Catalog Number: 65005

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid white
Molecular Weight	1171.18
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

EDTA succinimidyl ester is an excellent building block to develop site-specific ethylenediaminetetraacetic acid (EDTA) probes for variety of biomedical applications. It readily reacts with an amino compound under extremely mild conditions. EDTA is the most commonly used chelator for complexing transitional metal ions. EDTA chelation therapy uses the chelating agent, EDTA, to remove heavy metals and minerals from the body. It involves injecting EDTA into the bloodstream, where it binds to these substances, allowing them to be excreted through urine. EDTA is approved for treating lead poisoning, but it's also used in other conditions, including coronary artery disease and arthritis.