

EDTA Maleimide

Catalog Number: 65006

Unit Size: 5 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	24 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	690.66
------------------	--------

Soluble In	DMF
------------	-----

Spectral Properties

Excitation Wavelength	N/A
-----------------------	-----

Emission Wavelength	N/A
---------------------	-----

Applications

EDTA maleimide is an excellent building block to develop site-specific ethylenediaminetetraacetic acid (EDTA) probes for variety of biomedical applications. It readily reacts with a thiol 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.