

XFD594 C5 Maleimide

Catalog Number: 1891

Unit Size: 1 mg

Product Details

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

Expiration Date	12 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid purple
------------	--------------

Molecular Weight	1001.11
------------------	---------

Soluble In	DMSO
------------	------

Spectral Properties

Excitation Wavelength	590 nm
-----------------------	--------

Emission Wavelength	618 nm
---------------------	--------

Applications

N/A