

Phe-AMC

Catalog Number: 13423

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	322.36
------------------	--------

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

Spectral Properties

Excitation Wavelength	341 nm
-----------------------	--------

Emission Wavelength	441 nm
---------------------	--------

Applications

Phe-AMC is a protease inhibitor that inhibits the activity of serine and cysteine proteases. It binds to the active site of these enzymes and blocks their function. It is also used for monitoring the activity of α -chymotrypsin. It has also been shown to have biological functions such as the prevention of carcinogenesis.