

Cys-AMC

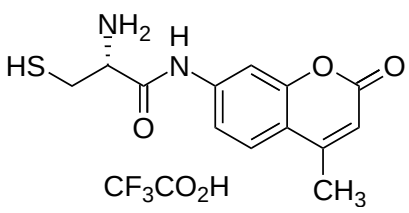
Catalog Number: 13493

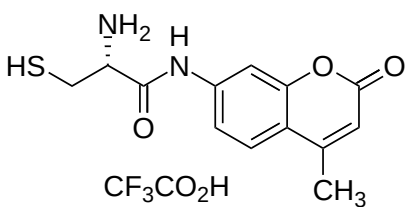
Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	N/A
Molecular Weight	392.35
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	341 nm
Emission Wavelength	441 nm

Applications

Cys-AMC is a sensitive fluorogenic protease substrate. The non-fluorescent substrate generates a bright blue fluorescent AMC product that has an emission spectra that can be easily detected with the DAPI filter set. It can be used for monitoring protease activity, making it a valuable research tool in enzymology, inhibitor screening, and mechanistic protease studies.