

Glu-AMC

Catalog Number: 13494

Unit Size: 5 mg

Product Details

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

Expiration Date 12 months upon receiving

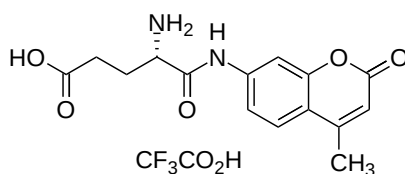
Chemical Properties

Appearance N/A

Molecular Weight 418.32

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength 341 nm

Emission Wavelength 441 nm

Applications

Glu-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.