

Ac-Leu-AMC

Catalog Number: 13444

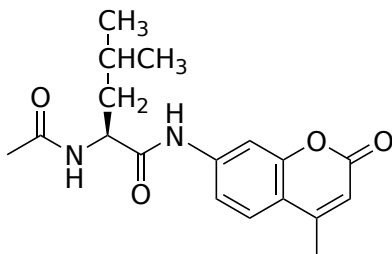
Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid off-white
Molecular Weight	330.38
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	341 nm
Emission Wavelength	441 nm

Applications

Ac-Leu-AMC is a fluorogenic substrate for leucine aminopeptidase. Upon the hydrolysis by proteases, Leu-AMC releases the strongly fluorescent AMC fluorophore that can be readily detected with Ex/Em = 350/430 nm. Leucine aminopeptidase is an extracellular enzyme found abundantly in natural aquatic systems. It is one of the various enzymes used by bacteria for protein hydrolysis. Ac-L-Leucine-7-amido-4-methylcoumarin might be used as a substrate in fluorescence-based soil assays to determine leucine aminopeptidase activity of both *Plasmodium falciparum* M1 (PfA-M1) and PfA-M17 enzymes.