

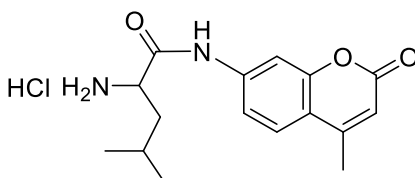
Product Information Sheet

Ordering Information

Product Number:	13445
Product Name:	Leu-AMC
Unit Size:	10 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 324.81



Chemical Structure:

Soluble in:	DMSO
Excitation Wavelength:	351 nm
Emission Wavelength:	430 nm

Application Notes

Leu-AMC is a fluorogenic substrate for leucine aminopeptidase. The weakly fluorescent substrate generates a bright blue fluorescent AMC product that has Ex/Em = 351/430 nm, and can be easily detected with the DAPI filter set.