

## D-VLK-AMC [D-Val-Leu-Lys-AMC]

Catalog number: 13201 Unit size: 5 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Off-white solid
Molecular Weight	743.69
Soluble In	DMSO
Chemical Structure	

Spectral Properties		
Excitation Wavelength	341 nm	
Emission Wavelength	441 nm	

## Applications

D-VLK-AMC is a sensitive, highly specific fluorescent substrate for plasmin. It can be used with a fluorescence microscope, fluorescence microplate or flow cytometer.