

AFC [7-Amino-4-trifluoromethylcoumarin] *CAS 53518-15-3*

Catalog number: 56, 57, 58 Unit size: 1 g, 5 g, 25 g

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	24 months upon receiving
Chemical Properties	
Appearance	Yellow solid
Molecular Weight	229.15
Soluble In	DMSO
Chemical Structure	H_2N H_2N G
Spectral Properties	
Excitation Wavelength	376 nm

Excitation Wavelength	3	76 nm
Emission Wavelength	48	82 nm

Applications

This compound is the precursor for a variety of AFC enzyme substrates. It is also used as a calibration standard for AFC-based enzyme substrates. Compared to AMC enzyme substrates, fluorogenic AFC substrates generate AFC that has longer absorption and fluorescence wavelength that AMC.