

AMP-Fluorescein conjugate calibrator

Catalog number: 13610 Unit size: 100 nmol

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Orange solid
Molecular Weight	~800
Soluble In	DMSO
Chemical Structure	$HO_{P}O_{OH}O_{OH}O_{O}$
Spectral Properties	
Excitation Wavelength	498 nm
Emission Wavelength	517 nm

Applications

AMP-Fluorescein conjugate calibrator is designed to use with our green fluorescent FAM-cAMP PDE IV substrate (#Cat: 13602), a cAMP derivative that is a specific substrate for phosphodiesterase (PDE) IV. This AMP-Fluorescein conjugate calibrator is the hydrolyzed product of FAM-cAMP PDE IV substrate by PDE IV. It is used to calibrate the hydrolysis rate of FAM-cAMP PDE IV substrate by PDE IV.