

## Product Information Sheet

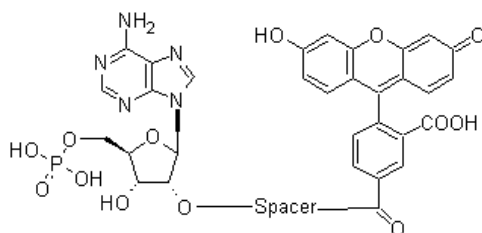
### Ordering Information

Product Number:	13610
Product Name:	AMP-Fluorescein conjugate calibrator
Unit Size:	100 nmol
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: ~800

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	492 nm
Emission Wavelength:	515 nm

### Application Notes

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 (#Cat: 13602) by PDE IV. It is used to calibrate the hydrolysis rate of FAM-cAMP PDE IV substrate (#Cat: 13602) by PDE IV.