

GMP-Fluorescein conjugate calibrator

Catalog number: 13611 Unit size: 100 nmol

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Solid

Molecular Weight ~800

Soluble In DMSO

Chemical Structure

Spectral Properties

Excitation Wavelength 498 nm

Emission Wavelength 517 nm

Applications

GMP-Fluorescein conjugate calibrator is designed to use with our green fluorescent FAM-cGMP PDE V substrate (#Cat: 13604), a cGMP derivative that is a specific substrate for phosphodiesterase (PDE) V. This GMP-Fluorescein conjugate calibrator is the hydrolyzed product of FAM-cGMP PDE V substrate by PDE V. It is used to calibrate the hydrolysis rate of FAM-cGMP PDE V substrate by PDE V.