

## Mca-YVADAPK(Dnp)-OH

Catalog number: 13557 Unit size: 1 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Solid
Molecular Weight	1145.15
Soluble In	DMSO
Spectral Properties	
Excitation Wavelength	322 nm
Emission Wavelength	381 nm

## Applications

Mca-YVADAPK(Dnp)-OH is a fluorogenic peptide substrate for detecting caspase-1/interleukin-converting enzyme (ICE) and angiotensin Iconverting enzyme-2 (ACE-2). Its N-terminal contains a highly fluorescent 7-methoxycoumarin (MCA) fluorophore that is efficiently quenched by non-fluorescent 2,4-dinitrophenyl group on lysine residue via fluorescence resonance energy transfer (dark FRET). It can be used to measure the activities of caspase-1/interleukin-converting enzyme (ICE) and angiotensin I-converting enzyme-2 (ACE-2) that are capable of cleaving an amide bond between the fluorescent group (MCA) and the quencher group (DNP), resulting an increase in fluorescence.