# Mca-APK(Dnp)

#### **ACE2** substrate

Fluorogenic substrate for angiotensin converting enzyme-2 (ACE2), a zinc-dependent carboxypeptidase with 42% amino acid sequence identity to ACE1. Unlike Mca-YVADAPK(Dnp), this substrate is not cleaved by caspases. Mca fluorescence is quenched by the Dnp group until cleavage (at Pro-Lys) separates them. Ex: 328 nm, Em: 393 nm, although the following Ex/Em have also been used: 320-340/400-420. This highly quenched (99.6%) substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic Mca control peptide (BML-P127).

Citations: 25

**View Online »** 

**Ordering Information** 

**Order Online** »

BML-P163-0001

1mg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Stock solutions in

DMSO stable at -20°C

**Handling** Protect from light and moisture.

Long Term Storage -20°C

**Shipping** Blue Ice

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name ACE2 substrate (fluorogenic)

**Appearance** Yellow powder.

Formula  $C_{32}H_{37}N_6O_{12}$ 

**Identity** Determined by MS.

**MW** 696.7

Purity ≥95% (HPLC)

Sequence Mca-Ala-Pro-Lys(Dnp) [Mca=(7-methoxycoumarin-4-

yl)acetyl;Dnp=2,4-dinitrophenyl]

Solubility Soluble in DMSO or 10% Ammonium Hydroxide (2 mg/ml).

**Source** Synthetic.

Last modified: May 29, 2024

