# Matrix metalloproteinase-9 (MMP-9) fluorometric drug discovery kit,

Fluorometric (also known as fluorimetric) Drug Discovery Kit, RED is a complete assay system designed to screen MMP-9 inhibitors using a quenched fluorogenic substrate OMNIMMP® RED: TQ3-GABA-Pro-Cha-Abu-Smc-His-Ala-Dab(6-TAMRA)-Ala-Lys-NH<sub>2</sub> [TQ3=quencher; GABA=4-aminobutyric acid; Cha=L-cyclohexylalanine; Abu=2-aminobutyric acid; Smc=S-methyl-L-cysteine; Dab=2,4-diaminobutyric acid; 6-TAMRA=6-tetramethylrhodamine]. TAMRA fluorescence is thoroughly quenched by the TQ3 group until cleavage by MMPs separates the two moieties.

The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of MMP-9, a potential therapeutic target. The compound NNGH is also included as a prototypic control inhibitor.

Matrix metalloproteinase-9 (MMP-9, gelatinase B, 92kDa type IV collagenase) is a member of the MMP family of extracellular proteases. These enzymes play a role in many normal and disease states by virtue of their broad substrate specificities. Targets of MMP-9 include native and denatured collagens, fibronectin, elastin, laminin, pro-TNF- $\alpha$ , and interleukins and their receptors. MMP-9 is secreted as a 92kDa proenzyme (as measured by SDS-PAGE), and activated by cleavage to forms 82kDa and smaller. MMP-9 is an important target for inhibitor screening due to its involvement in diseases such as alopecia, cancer, angiogenesis, and metastasis.

# **Ordering Information**

**Order Online** »

BML-AK306-0001

96 wells

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Gelatinase B, 92kDa Type IV collagenase

**Application** Activity assay, Fluorescent detection, HTS

**Contents** 1 vial MMP-9 enzyme

1 vial Substrate (OMNIMMP<sup>®</sup> RED) 1 vial 6'-TAMRA calibration standard

1 vial control inhibitor (NNGH)1 bottle (20 ml) assay buffer1 black 96-well microplate

Instructions

Technical Info / Product Notes NCBI Reference Sequence: NM 004994

The OMNIMMP<sup>®</sup> RED substrate offers key advantages over other MMP substrates.

- Emission at the red end of the spectrum (576 nm after excitation at 545 nm) avoids the interference at lower wavelengths often exhibited by screening compounds, and by substances commonly found in biological samples and tissue culture medium.
- 2. MMP substrate peptides display poor aqueous solubility, often with  $K_m$ s near their limits of solubility, making enzyme and inhibitor kinetics difficult. MMP  $K_m$ s for OMNIMMP<sup>®</sup> RED substrate are well below its solubility limit.
- 3. OMNIMMP<sup>®</sup> RED is avidly cleaved by MMPs, with *k* <sub>cat</sub>/K<sub>m</sub>s in the range of ~10<sup>4</sup>-10<sup>6</sup> M<sup>-1</sup>sec<sup>-1</sup>.
   4. The ultra-strong fluorescence of OMNIMMP<sup>®</sup> RED
- 4. The ultra-strong fluorescence of OMNIMMP® RED allows for substrate concentrations much lower than the  $K_m$ , a condition generally desirable in inhibitor screening assays.

UniProt ID P14780

Last modified: May 29, 2024



info-

eu@enzolifesciences.com