

# Matrix metalloproteinase (MMP) inhibitor profiling kit, fluorometric RED

The MMP inhibitor profiling kit, fluorometric (also known as fluorimetric) RED is a complete assay system designed to examine the specificity of inhibitors against a panel of ten matrix metalloproteinase enzymes, 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 compound NNGH is also included as a prototypic control inhibitor.

The matrix metalloproteinases, or MMPs, are extracellular proteases that function at a neutral pH to cleave a wide variety of substrates. These include basement membrane and extracellular matrix components, growth and death factors, cytokines, and cell and matrix adhesion molecules. The broad range of substrate specificities and expression patterns of MMPs results in their involvement in many different processes, both normal and pathological. Aberrant expression has been noted in cancer, angiogenesis, arthritis, inflammation, periodontal disease, emphysema, multiple sclerosis, pre-eclampsia, and chronic wounds, among others. The general structure of an MMP protein consists of a pre domain to direct secretion from the cell, a pro domain, a catalytic domain, and a C-terminal hemopexin domain. The catalytic site involves a coordinately-bound zinc ion. The inactive, or zymogen, form of the enzyme is activated by disruption of one of the coordinate bonds, usually via proteolytic removal of the pro domain.

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-AK308-0001

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

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

**Application Notes** Designed to examine the specificity of inhibitors against a panel of ten matrix metalloproteinase enzymes, using a quenched fluorogenic peptide.

**Contents**

- 10 vials (10 MMP enzymes)
- 1 vial Substrate (OMNIMMP® RED)
- 1 vial 6'-TAMRA calibration standard
- 1 vial control inhibitor (NNGH)
- 2 bottles (20 ml each) assay buffer
- 1 black 96-well microplate
- Instructions

**Technical Info / Product Notes** **NCBI Reference Sequence:** NM\_002429, NM\_004530, NM\_002422, NM\_002423, NM\_002424, NM\_004994, NM\_002425, NM\_002426, NM\_002427, NM\_004995

The OMNIMMP® RED substrate offers key advantages over other MMP substrates.

1. 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® RED substrate are below its solubility limit.
3. In addition to the efficient binding as exhibited by low  $K_m$ s, OMNIMMP® RED is avidly cleaved by MMPs, with  $k_{cat}/K_m$ s in the range of  $10^4$ - $10^6$  M<sup>-1</sup>sec<sup>-1</sup>.
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.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)