

# Matrix

## metalloproteinase-7 (MMP-7) fluorometric drug discovery kit,

### RED

The MMP-7 Fluorometric (also known as fluorimetric) Drug Discovery Kit, RED is a complete assay system designed to screen MMP-7 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-7, a potential therapeutic target. The compound NNGH is also included as a prototypic control inhibitor.

Matrix metalloproteinase-7 (MMP-7, matrilysin, pump-1) 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-7 include collagen, osteopontin, pro-TNF- $\alpha$ , E-cadherin,  $\beta$ 4 integrin, and Fas ligand. MMP-7 is secreted as a 28kDa proenzyme (as measured by SDS-PAGE), and activated by cleavage to 19kDa. MMP-7 is an important target for inhibitor screening due to its involvement in diseases such as cancer.

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	Matrilysin, Pump
Application	Activity assay, Fluorescent detection, HTS
Application Notes	Designed to screen MMP-7 inhibitors using a quenched fluorogenic peptide.
Contents	1 vial MMP-7 enzyme 1 vial substrate (OMNIMMP <sup>®</sup> RED) 1 vial 6'-TAMRA calibration standard 1 vial control inhibitor (NNGH) 1 bottle (20 ml) assay buffer 1 black 96-well microplate Instructions

The OMNIMMP<sup>®</sup> 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<sup>®</sup> RED substrate are below its solubility limit.
3. In addition to the efficient binding as exhibited by low  $K_m$ s, OMNIMMP<sup>®</sup> 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<sup>®</sup> RED allows for substrate concentrations much lower than the  $K_m$ , a condition generally desirable in inhibitor screening assays.

UniProt ID

P09237

Last modified: May 29, 2024



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-  
be@enzolifesciences.com](mailto:info-be@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)