

Matrix metalloproteinase-2 (MMP-2) fluorometric drug discovery kit,

RED

The MMP-2 Fluorometric (also known as fluorimetric) Drug Discovery Kit, RED is a complete assay system designed to screen MMP-2 inhibitors using a quenched fluorogenic substrate OMNIMMP® RED: TQ3-GABA-Pro-Cha-Abu-Smc-His-Ala-Dab(6-TAMRA)-Ala-Lys-NH₂ [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-2, a potential therapeutic target. The compound NNGH is also included as a prototypic control inhibitor.

Matrix metalloproteinase-2 (MMP-2, gelatinase A, 72kDa 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-2 include native and denatured collagens, fibronectin, elastin, laminin-5, pro-TNF- α , and neurocan. MMP-2 is secreted as a 72kDa proenzyme (as measured by SDS-PAGE), and activated by cleavage to 62 and 59kDa. MMP-2 is an important target for inhibitor screening due to its involvement in diseases such as atherosclerosis, and cancer growth, angiogenesis, and metastasis.

Ordering Information

[Order Online »](#)

BML-AK302-0001	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 A, 72 kDa Type IV collagenase
Application	Activity assay, Fluorescent detection, HTS
Application Notes	Designed to screen MMP-2 inhibitors using a quenched fluorogenic peptide.
Contents	1 vial MMP-2 enzyme 1 vial substrate (OMNIMMP® 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[®] 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.

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⁻¹ sec⁻¹.

3. 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

P08253

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)