

Matrix

Metalloproteinase-14 (MMP-14) fluorometric drug discovery kit

A QUANTIZYME® Assay System. The MMP-14 Fluorometric (also known as fluorimetric) Drug Discovery Kit is a complete assay system designed to screen for inhibitors of matrix metalloproteinase-14 (MMP-14, membrane-type matrix metalloproteinase 1, MT1-MMP) using a quenched fluorogenic peptide: OMNIMMP® fluorogenic substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH. Mca fluorescence is quenched by the Dpa group until cleavage by MMPs at the Gly-Leu bond separates the two moieties. The assays are performed in a convenient 96-well microplate format. Also included are active enzyme, a prototypic control inhibitor (NNGH³), and a detailed instruction booklet.

Citations: 1

[View Online »](#)

Ordering Information

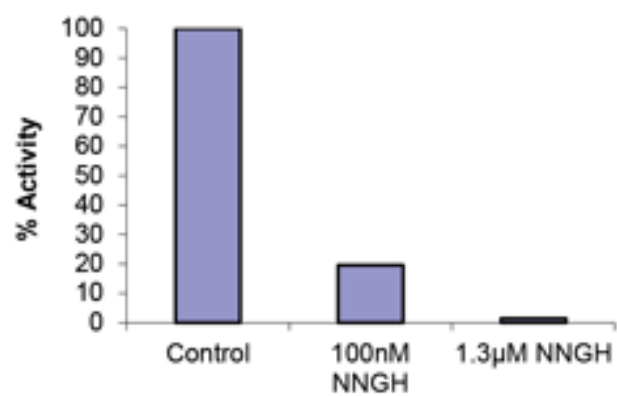
[Order Online »](#)

BML-AK417-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Inhibition of MMP-14 by NNGH



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Membrane type matrix metalloproteinase 1, MT-1 MMP

Application Activity assay, Fluorescent detection, HTS

Contents MMP-14 enzyme (human) (recombinant), substrate, calibration standard, Assay Buffer, 96-well white microplate.

UniProt ID P45452

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)