

Matrix

Metalloproteinase-8

(MMP-8) fluorometric

drug discovery kit

A QUANTIZYME® Assay System. The MMP-8 Fluorometric (also known as fluorimetric) Drug Discovery Kit is a complete assay system designed to screen inhibitors of matrix metalloproteinase-8 (MMP-8, neutrophil collagenase, collagenase-2) using a quenched fluorogenic peptide: OMNIMMP® fluorogenic substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH₂. 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.

Ordering Information

[Order Online »](#)

| | |
|----------------|----------|
| BML-AK415-0001 | 96 wells |
|----------------|----------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Neutrophil collagenase, Collagenase-2

Application Activity assay, Fluorescent detection, HTS

Contents MMP-8 enzyme (human) (recombinant), substrate, calibration standard, inhibitor, assay buffer, 96-well white microplate.

UniProt ID P22894



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)