

Matrix

Metalloproteinase-7 (MMP-7) colorimetric drug discovery kit

A QUANTIZYME® Assay System. The MMP-7 Colorimetric Drug Discovery Kit is a complete assay system designed to screen inhibitors of matrix metalloproteinase-7 (MMP-7, matrilysin) using a thiopeptide as a chromogenic substrate. The assays are performed in a convenient 96-well microplate format. Included are active enzyme, assay buffer, a prototypic control inhibitor (NNGH), and a detailed instruction booklet.

Ordering Information

[Order Online »](#)

BML-AK406-0001	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

Matrilysin, Pump

Application

Activity assay, Colorimetric detection, HTS

Contents

MMP-7 enzyme (human) (recombinant), substrate, inhibitor, assay buffer, 96-well microplate.

UniProt ID

P09237

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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