

# MMP Inhibitor profiling kit, colorimetric

A QUANTIZYME® Assay System Complete assay system designed to examine the specificity of inhibitors against a panel of ten matrix metalloproteinase enzymes. The assays are performed in a convenient 96-well microplate format using a chromogenic substrate<sup>1,2</sup> detected at 412nm. Included are active MMP enzymes, assay buffer, a prototypic control inhibitor (NNGH<sup>3</sup>), and a detailed instruction booklet.

## Ordering Information

[Order Online »](#)

|                       |          |
|-----------------------|----------|
| <b>BML-AK015-0001</b> | 96 wells |
|-----------------------|----------|

## Manuals, SDS & CofA

[View Online »](#)

- Complete assay system
- Optimized to detect the activity of a variety of highly active MMP enzymes
- Colorimetric detection, High throughput screening
- Convenient 96 well format

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Activity assay, Colorimetric detection, HTS

**Contents** MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-10, MMP-12, MMP-13, MMP-14 enzyme (human) (recombinant), substrate, inhibitor, assay buffer, 96-well microplate.

**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)