

Matrix

Metalloproteinase-2 (MMP-2) fluorometric drug discovery kit

A QUANTIZYME[®] Assay System. The MMP-2 Fluorometric (also known as fluorimetric) Drug Discovery Kit is a complete assay system designed to screen inhibitors of matrix metalloproteinase-2 (MMP-2, gelatinase A) 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. The kit is useful to screen inhibitors of MMP-2, a potential therapeutic target. Included are active enzyme, assay buffer, a prototypic control inhibitor (NNGH³), and a detailed instruction booklet.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK409-0001

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	Gelatinase A, 72 kDa Type IV collagenase
Application	Activity assay, Fluorescent detection, HTS
Contents	MMP-2 enzyme (human) (recombinant), substrate, calibration standard, inhibitor, assay buffer, 96-well white microplate.
UniProt ID	P08253

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