

MMP substrate (chromogenic)

Sensitive chromogenic MMP substrate

Chromogenic substrate for continuous spectrophotometric assay of most matrix metalloproteinases, and TACE (ADAM17). The MMP cleavage site peptide bond is replaced by a thioester bond in this peptide. Hydrolysis of this bond by an MMP produces a sulfhydryl group, which reacts with DTNB [5,5'-dithiobis(2-nitrobenzoic acid), Ellman's Reagent] to form 2-nitro-5-thiobenzoic acid, which can be detected by its absorbance at 412 nm.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P125-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After reconstitution, store at -20°C or -70°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Matrix metalloproteinase substrate, Ac-Pro-Leu-Gly-[2-mercapto-4-methyl-pentanoyl]-Leu-Gly-OC ₂ H ₅
Application Notes	Can be used for assay with human collagen cells.
Formulation	Lyophilized.
MW	655.9
Purity	≥97% (HPLC)
Sequence	Ac-Pro-Leu-Gly-[2-mercapto-4-methyl-pentanoyl]-Leu-Gly-OC ₂ H ₅
Solubility	Soluble in DMF (100mM) or DMSO (100mM).



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