

MMP substrate (fluorogenic)

Fluorogenic substrate for several matrix metalloproteinases (MMPs). It is cleaved by MMP-1, MMP-3, MMP-7, MMP-8, MMP-9, MMP-11, MMP-12, MMP-13, and ADAM9/MDC9, but only slightly by MMP-2 and not at all by ADAM17/TACE. Ex/Em= 340nm/440nm, although 365/450 is also appropriate. Highly quenched, with strong fluorescence by Nma at 440nm once cleavage separates it from the aromatic Dnp moiety. Compatible with most fluorometers.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P128-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Matrix metalloproteinase substrate, Dnp-Pro- β -cyclohexyl-Ala-Gly-Cys(Me)-His-Ala-Lys(Nma)-NH ₂
Appearance	lyophilized solid
MW	1077.4
Purity	≥98% (HPLC)
Sequence	Dnp-P(Cha)G(CMe)HA(KNma)-amide
Solubility	Soluble in DMSO (at least 20mM).



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)