

MMP substrate α (fluorogenic)

Fluorogenic substrate for several matrix metalloproteinases (MMPs). It is cleaved by MMP-1 ($k_{\text{cat}}/K_m = 3.3 \times 10^4 \text{ M}^{-1} \text{ s}^{-1}$), MMP-2 ($k_{\text{cat}}/K_m = 5.8 \times 10^4$), MMP-3 ($k_{\text{cat}}/K_m = 2.2 \times 10^4$), MMP-7 ($k_{\text{cat}}/K_m = 1.2 \times 10^4$), MMP-8, MMP-9 ($k_{\text{cat}}/K_m = 6.9 \times 10^4$), and MMP-26 ($k_{\text{cat}}/K_m = 4.9 \times 10^4$) Ex/Em=280nm/346nm, although 283/350,360 are also appropriate. Fluorescence at 346 nm by tryptophan occurs once cleavage separates it from the quenching aromatic Dnp moiety. MMPs cleave this peptide between Gly-Leu.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P131-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Matrix metalloproteinase substrate α
CAS	121282-17-5
Formula	$C_{45}H_{64}N_{14}O_{11}$
MW	977.09
Purity	≥94% (HPLC)
Sequence	Dnp-Pro-Leu-Gly-Leu-Trp-Ala-D-Arg-NH ₂ (Dnp=2,4-dinitrophenyl)

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-
fr@enzolifesciences.com](mailto:info-fr@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)