

Dnp-Pro-Leu-Gly-Met-Trp-Ser-Arg-OH

Substrate for MMP-2 and -9

Fluorogenic substrate for MMP-2 and MMP-9. Excitation: 280nm; emission 360nm.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-120-M005	5mg
------------------	-----

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow solid

Formula $C_{44}H_{61}N_{13}O_{13}S$

MW 1012.1

Purity ≥95% (HPLC)

Sequence 2,4-Dinitrophenyl-Pro-Leu-Gly-Met-Trp-Ser-Arg-OH

Solubility Soluble in DMSO; dilute with distilled water or buffer.

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)