

Ac-Ala-Pro-Nle-Asp-OH

Highly specific inhibitor of the caspase-like peptidic activity of the 20S proteasome.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AW9485-0100

100µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Once dissolved use solutions immediately and do not store.
Handling	Do not add buffers containing free amino groups as these will react with the aldehyde moiety thus inactivating the compound.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Glassy lyophilized solid.
Formula	$C_{20}H_{32}N_4O_8$
MW	456.5
Purity	≥95% (Reverse-Phase HPLC)
Sequence	Ac-Ala-Pro-Nle-Asp-OH
Solubility	Soluble in water:acetonitrile containing 0.1% TFA (3:1) at a concentration of 1mg/mL.



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)