

Z-Pro-Nle-Asp-CHO

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

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|--------------------|----------------------------|
| Appearance | White to off-white solid. |
| Formulation | Lyophilized. |
| MW | 478.5 |
| Purity | ≥95% (HPLC) |
| Sequence | Z-Pro-Nle-Asp-H |
| Solubility | Soluble in water (1mg/ml). |

Last modified: May 29, 2024



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