

Gelonin

Type I ribosome inactivating protein (RIP). Depurinates RNA in ribosomes, thus inhibiting protein synthesis in eukaryotic cells, which results in cell death. Widely used to construct immunotoxins composed of cell-targeted antibodies. As a type I RIP it lacks the lectin subunit and is practically non-toxic to intact cells. Inhibits cell-free protein synthesis in reticulocyte assay with globin messenger ($IC_{50}=4.5ng/ml$).

Citations: 26

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-350-150-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

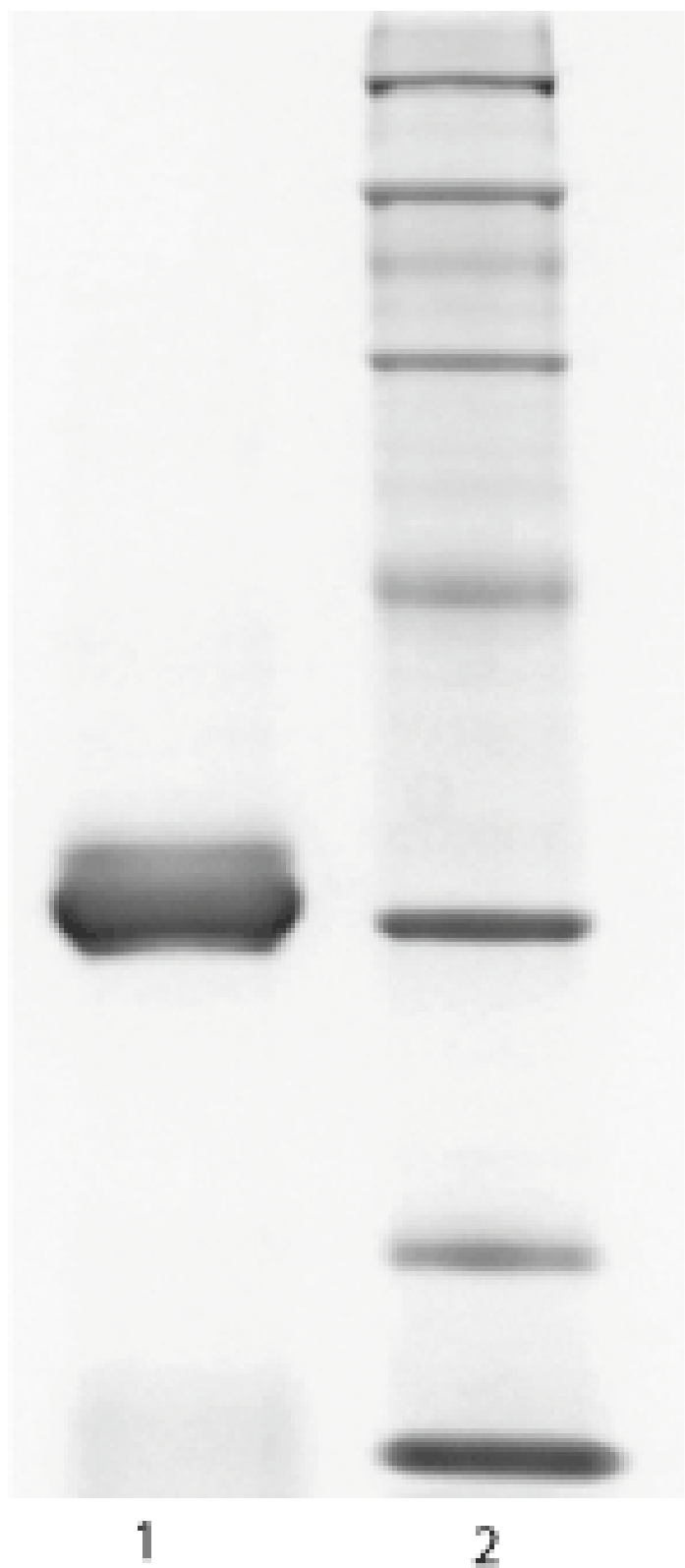


Figure: SDS-PAGE (12% polyacrylamidgel)

Lane 1: Gelonin, 5 μ g

Lane 2: Proteinmarker (200-6,5 kDa)

Handling & Storage

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

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

CAS 75037-46-6

Concentration ~2mg/ml

Formulation Liquid. In 5mM phosphate buffer, pH 6.5, containing 5mM EDTA and 120mM sodium chloride.

MW ~30kDa

Purity ≥95%

Source Isolated from *Gelonium multiflorum*.

UniProt ID P33186

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-
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)