

RGD peptide

Integrin ligand

Induces apoptosis in various cell types and cell lines at 1mM by a mechanism involving cellular uptake and possible direct stimulation of procaspase-3 autoprocessing. Does not induce apoptosis in the breast cancer line MCF-7, in which there is a functional deletion of the procaspase-3 gene. Contains the motif RGD (Arg-Gly-Asp) which is found in many integrin ligands. Inhibits cell adhesion of COS-7 cells to fibronectin 80-90% at 1mM.

Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P700-0005	5mg
---------------	-----

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 GRGDNP

Appearance Solid.

CAS 114681-65-1

Formulation Lyophilized.

MW 614.6

Purity ≥97% (HPLC)

Sequence Gly-Arg-Gly-Asp-Asn-Pro

Solubility Soluble in water (>50mg/ml).

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)