

Exendin-4

GLP-1 receptor agonist

Exendin-4 is a 39 aa polypeptide and incretin mimetic which was first isolated from Gila monster saliva. The synthetic form, Exenatide, is a glucagon like peptide-1 (GLP-1) agonist approved for the treatment of diabetes mellitus type II. Research shows Exendin-4 is involved in glucose-dependent enhancement of insulin secretion, reduction of hyperglycemia, glucose-dependent suppression of inappropriately high glucagon secretion, slowing of gastric emptying, and reduction of food intake, often with body weight reduction or blunting of weight gain.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT111-0001	1mg
ENZ-PRT111-0500	500µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Exenatide

Appearance White to off-white solid.

CAS 141758-74-9

Formulation Lyophilized.

MW 4186.57

Purity ≥98% (HPLC)

Sequence His-Gly-Glu-Gly-Thr-Phe-Thr-Ser-Asp-Leu-Ser-Lys-Gln-Met-Glu-Glu-Glu-Ala-Val-Arg-Leu-Phe-Ile-Glu-Trp-Leu-Lys-Asn-Gly-Gly-Pro-Ser-Ser-Gly-Ala-Pro-Pro-²Phe-Ser-NH₂

Solubility Soluble in water or DMSO.

Source Synthetic.



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)