

# Glucagon like peptide 1 (human) (7-36) amide

A predominant cleaved form of pro-glucagon expressed in the intestine. GLP-1 [7-36]amide is promptly released from the gut after oral ingestion of a meal where it slows gastric emptying and exerts a powerful insulinotropic action leading to elevated serum insulin levels. GLP-1 [7-36] amide promotes satiety and reduces food intake in patients with diabetes mellitus type 2 and has been suggested to have therapeutic potential in the treatment of NIDDM. It has been reported that human pancreatic endocrine tumors contain molecular forms of GLP-1 which resemble those expressed in the small intestine, rather than in their tissue of origin.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-GP1438-0100	100µg
-----------------	-------

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	GLP 1 (7-36) amide
Appearance	Lyophilized solid.
MW	3297.7
Sequence	H-His-Ala-Glu-Gly-Thr-Phe-Thr-Ser-Asp-Val-Ser-Ser-Tyr-Leu-Glu-Gly-Gln-Ala-Ala-Lys-Glu-Phe-Ile-Ala-Trp-Leu-Val-Lys-Gly-Arg-NH <sub>2</sub>
Solubility	Soluble in water at 1mg/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-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)