

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

A predominant cleaved form of pro-glucagon found in the pancreas and central nervous system. It has been reported that human pancreatic endocrine tumours contain molecular forms of GLP-1 which resemble those expressed in the small intestine, rather than in their tissue of origin. Within the CNS, GLP-1 powerfully inhibits feeding in fasting rats, an action which is blocked by prior administration of the GLP-1 receptor antagonist exendin [9-39]. This observation led to the suggestion that central GLP-1 is a physiological mediator of satiety and has projected GLP-1 into the obesity control debate. Glucagon-like peptide (GLP) 1 [1-36] also serves as a control peptide for Anti-Glucagon-like Peptide (GLP) 1 (BML-GA1176). Blocks immunoreactivity at 10 nmol per ml diluted antibody.

**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 (1-36) amide, Blocking peptide for Glucagon like peptide 1 pAb (Prod. No. BML-GA1176)
Appearance	Lyophilized solid.
MW	4111.5
Purity	≥95%
Sequence	H-His-Asp-Glu-Phe-Glu-Arg-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-amide
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)