

# Exendin-4 monoclonal antibody (20)

Exendin-4 is a 39-amino-acid peptide found in venom from the Gila monster *Heloderma suspectum*. It is a member of the glucagon-secretin family of peptide hormones and neuropeptides. Exendin-4 is a potent agonist of the GLP-1 receptor, and hence a potent stimulator of insulin secretion.

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

BPD-ABS-012-20-1	1mg
BPD-ABS-012-20-04	400µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application ELISA

Application Notes Detects a band of ~4kDa by Western blot.

Clone 20

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic exendin-4 from Gila monster *Heloderma suspectum*.

Isotype IgG1κ

Purity Detail Protein A/G-affinity purified.

Species Reactivity Gila monster (*Heloderma suspectum*)

Specificity Recognizes Gila monster *Heloderma suspectum* exendin-4 which has a high degree of homology with GLP-1 found in mammals.

UniProt ID P26349

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)