

Exendin-4 monoclonal antibody (35) (biotin conjugate)

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](#).

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 35

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14mM sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic exendin-4.

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](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com