

# Exendin-4 monoclonal antibody (35)

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-012-35-04	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 35

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](#)

.