

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 <i>Heloderma suspectum</i> .
Isotype	IgG1k
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Gila monster (<i>Heloderma suspectum</i>)
Specificity	Recognizes Gila monster <i>Heloderma suspectum</i> 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