

Ghrelin monoclonal antibody (121) (biotin conjugate)

Ghrelin is an orexigenic (food intake stimulating) 27 or 28 amino-acid gut-brain peptide hormone produced by endocrine cells in the stomach, pancreatic islets and certain hypothalamic neurons. Two forms are produced by alternative mRNA splicing; Gln13 or 14 is missing from the second form. "Active" ghrelin is acylated (octanoylated) at Ser3; this form stimulates growth hormone secretion via GHS receptors and increases appetite. The des-acyl form does not have these effects but may have other actions. Plasma responses of total ghrelin to food may reflect those of active ghrelin.

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

Ordering Information

[Order Online »](#)

| | |
|----------------------|-------|
| BPD-ABS-052-121B-015 | 150µg |
|----------------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Clone 121

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

GenBank ID 51738

Host Mouse

Immunogen Synthetic peptide corresponding to aa 14-28 of human ghrelin.

Isotype IgG2bk

Purity Detail Protein A purified.

Species Reactivity Cat, Dog, Human, Mouse, Rat

UniProt ID Q9UBU3

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)