

Ghrelin monoclonal antibody (121)

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](#).

Citations: 1

[View Online »](#)

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, WB
Clone	121
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M 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.
Source	From culture supernatant.
Species Reactivity	Cat, Dog, Human, Mouse, Rat
UniProt ID	Q9UBU3
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