Oxyntomodulin monoclonal antibody (46)

Oxyntomodulin is a peptide hormone of 37 amino-acid residues (29-amino acid sequence of glucagon, with an 8-amino acid carboxy-terminal extension) produced by the L cells of the gut. It modulates gastric acid secretion from the oxyntic cells of the stomach, inhibits appetite and has glucose lowering effects. Oxyntomodulin is produced by cell-type-specific processing of proglucagon that forms as well glicentin, GLP-1 and GLP-2 in the L cells of the intestinal mucosa.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-064-46-1	1mg
BPD-ABS-064-46-04	400μg

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 46

Crossreactivity Does not cross-react with glucagon.

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

GenBank ID 2641

Host Mouse

Immunogen Synthetic oxyntomodulin 31-37 coupled to carrier.

Isotype IgG1ĸ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human, Mouse, Rat

Specificity Specific for C-terminus of human oxyntomodulin.

UniProt ID P01275

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com