GLP-1 (7-37, 7-36, amide, free NT) monoclonal antibody (10) (biotin conjugate)

Glucagon-like peptide 1 (7-36) amide (GLP-1 (7-36) amide) is the principal active form of GLP-1, the other being GLP-1 (7-37). GLP-1 is a peptide hormone of the glucagon family, produced by the L cells of the intestinal mucosa from the same prohormone as glucagon. The active forms are potent stimulators of glucose-dependent insulin secretion. The sequence of GLP-1 is fully conserved in all mammalian species examined so far.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

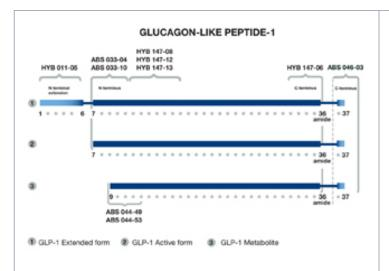
Order Online »

BPD-ABS-033-10B-015

150µg

Manuals, SDS & CofA

View Online »



Schematic glucagon-like peptide-1 (GLP-1) showing active, extended and metabolite forms along with select GLP-1 monoclonal antibodies.

Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Glucagon like peptide 1

Application ELISA

Clone 10

Crossreactivity Shows <0.2% cross-reactivity with GLP-1(9-36)amide,

glucagon and human GIP. Cross-reacts 0.25% with GLP-

1(1-37), approximately 6% with human GLP-2 and

approximately 50% with exendin-4.

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

0.14mM sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 7-17 of GLP-1.

lsotype lgG2aк

Purity Detail Protein A/G-affinity purified.

Species Reactivity Mammalian

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

France
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