

Glucagon like peptide 2 (human)

Glucagon-like peptide 2 is a 33- or 34-amino acid peptide located at the C-terminus of, and cleaved from, *pro*-glucagon by prohormone convertase 3 (PC3). GLP-2 is a potent intestinotropic growth factor, the action of which is transduced by G protein-coupled membrane-bound receptors (GLP2-R) found in the gut. GLP-2Rs are also found in the rodent CNS. The main actions of GLP-2 appear to be crypt cell proliferation, inhibition of enterocyte apoptosis, enhancement of nutrient adsorption and certain reparative effects. Applications: antiserum adsorption control; standard or tracer for immunoassay.

Ordering Information

[Order Online »](#)

BML-GP1439-0100	100µg
------------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GLP 2

Appearance Lyophilized solid.

CAS 99120-49-7

MW 3922.4

Purity ≥90%

Sequence H-His-Ala-Asp-Gly-Ser-Phe-Ser-Asp-Glu-Met-Asn-Thr-Ile-Leu-Asp-Asn-Leu-Ala-Ala-Arg-Asp-Phe-Ile-Asn-Trp-Leu-Ile-Gln-Thr-Lys-Ile-Thr-Asp-Arg-OH

Solubility Soluble in dilute acetic acid at 1mg/mL.

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-fr@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