

Product Data Sheet

Product Name: Glucagon-like Peptide-2, GLP-2 (146-178), human

Catalog Number: AS-62070 (1 mg) Lot Number: See label on vial

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-

OH (3-letter code)

HADGSFSDEMNTILDNLAARDFINWLIQTKITD (1-letter code)

Molecular Weight: 3766.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Glucagon-like Peptide-2 is freely soluble in DMSO.

Storage: Glucagon-like Peptide-2 is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at – 20°C or lower.

Description: This is a fragment of human intestinal growth factor glucagon-like peptide 2, GLP2, containing amino acids 146 to 178. It is an intestinotrophic growth hormone that promotes many aspects of intestinal function, including enhancement of mucosal growth and promotion of nutrient absorption. GLP-2 is a hormone that can rapidly improve intestinal epithelial barrier function. Ref: Cameron, H. et al. *JPET* 314, 214 (2005).

Related Products:

Name Cat # Size
GLP-2 (1-34), Glucagon-Like Peptide-2 (1-34), human AS-62068 1 mg
HADGSFSDEMNTILDNLAARDFINWLIQTKITDR

For Research Use Only