



Product Data Sheet

Product Name:	Glucagon-Like Peptide 1, GLP-1 amide, human	
Catalog Number:	AS-22460 (0.5 mg)	Lot Number: See label on vial
Sequence:	H-His-Asp-Glu-Phe-Glu-Arg-His-Ala-Glu-Gly-Thr-Phe-Thr-Ser-Asp-Val-Ser-Ser-Tyr-Leu-Glu-Gly-Gln-Ala-Ala-Lys-Glu-Phe-Ile-Ala-Trp-Leu-Val-Lys-Gly-Arg-NH ₂ (3-letter code) HDEFERHAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH ₂ (1-letter code)	
Molecular Weight:	4111.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Initially try to dissolve a small amount of Glucagon-Like peptide in water. If peptide does not dissolve in water, then use 1% NH₄OH.

Storage: Glucagon-Like Peptide 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 peptide, formed by cleavage of two basic residues, is an endogenous peptide secreted into the circulation by intestinal L cells. A potent glucose-dependent insulinotropic hormone, it has important actions on gastric motility, on the suppression of plasma glucagon levels and possibly on the promotion of satiety and stimulation of glucose disposal in peripheral tissues independent of the actions of insulin. It potentiates insulin secretion by pancreatic A-cells in a glucose-dependent manner and inhibits gastric emptying and acid secretion. Ref: Drucker, D. et al. *Proc. Natl. Acad. Sci. USA* **84**, 3434 (1987); Kieffer, T. and J. Habener, *Endo. Rev.* **20**, 876 (1999).

Related Products:

Name	Cat #	Size
Glucagon-Like Peptide 1, GLP-1 amide, human, FAM-labeled	AS-23606	0.5 mg
FAM-HDEFERHAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH ₂	AS-23607	1 mg
Biotin-Glucagon-Like Peptide 1, GLP-1 (7-36), amide, human	AS-23586	0.5 mg
Biotin-HAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH ₂		
Glucagon-Like Peptide 1, GLP-1 (7-36)-Lys(Biotin), amide, human	AS-23642	0.5 mg
HAEGTFTSDVSSYLEGQAAKEFIAWLVKGRK(Biotin)-NH ₂	AS-23643	1 mg

Glucagon-Like Peptide 1, GLP-1 (7-37)	AS-20761	0.5 mg
HAEGTFTSDVSSYLEGQAAKEFIAWLVKGRG	AS-20762	1 mg
Glucagon-Like Peptide 1, GLP-1 (7-36), amide, human	AS-22462	0.5 mg
HAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH ₂	AS-22463	1 mg
[FAM-Trp31]-GLP-1 (7-36)	AS-64488	0.5 mg
HAEGTFTSDVSSYLEGQAAKEFIA (Trp-S-FAM)		

For Research Use Only