

Product Data Sheet

Product Name: Glucagon-Like Peptide 1, GLP-1 (7-36), amide, human

Catalog Number: AS-22463 (1 mg) Lot Number: See label on vial

Sequence: H-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-NH2

(3-letter code)

HAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH2 (1-letter code)

Molecular Weight: 3298.7

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Glucagon-Like peptide is freely soluble in H₂O.

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 shares sequence with preproglucagon (78-107), amide, human.

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Human GLP-1 (7–36) amide (AnaSpec, San Jose, CA, USA) was infused at a constant rate of 1.5 pmol kg⁻¹ min⁻¹ via a mini-osmotic pump that delivers the solution continuously for up to 28 days (ALZET, model 2004; DURECT, Cupertino, CA, USA) as described previously. To initiate immediate pumping, we placed the pre-filled pumps in 0.8% saline for at least 40 h at 37°C before implantation. According to the manufacturer's guideline, GLP-1 was dissolved in 5% acetic acid buffer-Zhang, J. et al. *Diabetologia* 50, 1900 (2007). Published Citations:

Zhang, J. et al. *Diabetologia* 50, 1900 (2007).

Related Products:

Name	Cat #	Size
Biotin-Glucagon-Like Peptide 1, GLP-1 (7-36), amide, human	AS-23586	0.5 mg
(Biotin-HAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-NH2)		
Glucagon-Like Peptide 1, GLP-1 (7-36)-Lys(Biotin), amide, human	AS-23642	0.5 mg
	AS-23643	· ·
(HAEGTFTSDVSSYLEGQAAKEFIAWLVKGRK(Biotin)-NH2)	A3-23043	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 amide, human, FAM-labeled	AS-23606	0.5 mg
(FAM-HDEFERHAEGTFTSDVSSYLEGOAAKEFIAWLVKGR)	AS-23607	1 mg

For Research Use Only