



Product Data Sheet

Product Name:	Glucagon (1-29), bovine, human, porcine	
Catalog Number:	AS-22456 (0.5 mg) AS-22457 (1 mg)	Lot Number: See label on vial
Sequence:	H-His-Ser-Gln-Gly-Thr-Phe-Thr-Ser-Asp-Tyr-Ser-Lys-Tyr-Leu-Asp-Ser-Arg-Arg-Ala-Gln-Asp-Phe-Val-Gln-Trp-Leu-Met-Asn-Thr-OH (3-letter code) HSQGTFTSDYSKYLDSRRAQDFVQWLMNT (1-letter code)	
Molecular Weight:	3482.8	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Glucagon (1-29) peptide is freely soluble in water or 1% NH₄OH.

Storage: Glucagon (1-29) peptide 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.

Related Products:

Name	Cat #	Size
Biotin-Glucagon (1-29), bovine, human, porcine	AS-60274-05	0.5 mg
Biotin-HSQGTFTSDYSKYLDSRRAQDFVQWLMNT	AS-60274-1	1 mg
Biotin-Glucagon (1-29), bovine, human, porcine	AS-60273-1	1 mg
Biotin-HSQGTFTSDYSKYLDSRRAQDFVQWLMNT		

For Research Use Only