



Product Data Sheet

Product Name: Biotin-Gastrin (1-17)
Catalog Number: AS-60329-1 (1 mg) **Lot Number:** See label on vial
Sequence: Biotin-Glu-Gly-Pro-Trp-Leu-Glu-Glu-Glu-Glu-Glu-Ala-Tyr-Gly-Trp-Met-Asp-Phe-NH₂ (3-letter code)
Biotin-EGPWLEEEEEAYGWMDf-NH₂ (1-letter code)
Molecular Weight: 2343.6
% Peak Area by HPLC: ≥ 95
Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 100 µl 1% NH₄OH to 1 mg Biotin-Gastrin peptide. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20°C.

Storage: Biotin-Gastrin 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.

Related Products:

Name	Cat #	Size
Gastrin Releasing Peptide, human (VPLPAGGGTVLTKMYPRGNHWAVGHLM-NH ₂)	AS-24213 AS-24214	0.5 mg 1 mg
Big Gastrin-1, human (Pyr-LGPQGPPHLVADPSKKQGPWLEEEEEAYGWMDf-NH ₂)	AS-20747	0.5 mg

For Research Use Only