



Product Data Sheet

| | | |
|--------------------------------|---|-------------------------------|
| Product Name: | GIP (1-42), human | |
| Catalog Number: | AS-61226-05 (5 mg) | Lot Number: See label on vial |
| Sequence: | H-Tyr-Ala-Glu-Gly-Thr-Phe-Ile-Ser-Asp-Tyr-Ser-Ile-Ala-Met-Asp-Lys-Ile-His-Gln-Gln-Asp-Phe-Val-Asn-Trp-Leu-Leu-Ala-Gln-Lys-Gly-Lys-Lys-Asn-Asp-Trp-Lys-His-Asn-Ile-Thr-Gln-OH (3-letter code) YAEGTFISDYSIAMDKIHQQDFVNWLLAQKGKKNWVKHNITQ (1-letter code) | |
| Molecular Weight: | 4983.6 | |
| Peptide Purity: | >95% | |
| Appearance: | Lyophilized white powder | |
| Peptide Reconstitution: | Using DMSO, reconstitute by adding 50-70 µl to 0.5 mg GIP. | |
| Storage: | GIP 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. | |

For Research Use Only