



Product Data Sheet

| | | |
|--------------------------|---|-------------------------------|
| Product Name: | Galanin, human | |
| Catalog Number: | AS-22431 (1 mg) | Lot Number: See label on vial |
| Sequence: | H-Gly-Trp-Thr-Leu-Asn-Ser-Ala-Gly-Tyr-Leu-Leu-Gly-Pro-His-Ala-Val-Gly-Asn-His-Arg-Ser-Phe-Ser-Asp-Lys-Asn-Gly-Leu-Thr-Ser-OH (3-letter code) GWTLNSAGYLLGPHAVGNHRFSFSDKNGLTS(1-letter code) | |
| Molecular Weight: | 3157.5 | |
| Peptide Purity: | >95% | |
| Appearance: | Lyophilized white powder | |

Peptide Reconstitution: Galanin peptide is freely soluble in water.

Storage: Galanin 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: Galanin, a 29 or 30 (in human) amino acid neuropeptide, is found in the central and peripheral nervous systems and displays several important physiological activities. These actions are mediated through distinct G protein-coupled receptors.

Published Citations:

- Ahn, H. et al. *Mol. Pharmacol.* **51**, 350 (1997).
- Mazarati, Andréy. et al. *J. Pharmacol.* **318**, 700 (2006).
- Misawa, K. et al. *Clin Cancer Res* **14**, 7604 (2008).
- Kanazawa, T. et al. *Clin Cancer Res* **15** 2222 (2009).

For Research Use Only