



Product Data Sheet

Product Name: Bombesin

Catalog Number: AS-20665 (1 mg) **Lot Number:** See label on vial

Sequence: Pyr-Gln-Arg-Leu-Gly-Asn-Gln-Trp-Ala-Val-Gly-His-Leu-Met-NH2
(3-letter code)
Pyr-QRLGNQWAVGHLM-NH2 (1-letter code)

Molecular Weight: 1620.9

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Bombesin peptide is freely soluble in water.

Storage: Bombesin 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 14-amino-acid bombesin peptide is homologous to the C terminus of Gastrin-Releasing Protein (GRP). It binds with high affinity to the GRP receptor, a member of the G-protein-coupled receptor family. Ref: Dietrich, JB. *Cell. Mol. Biol.* **40**, 731 (1994).

Published Citations:

Bogunovic, Milena. et al. *Am J Physiol Gastrointest Liver Physiol* **292**, G1770 (2007).
O'Donoghue, A. et al. *J Biol Chem* **283**, 29186 (2008).

Related Products:

Name	Cat #	Size
[Lys3]-Bombesin Pyr-QKLGNGWAVGHLM-NH2	AS-22838	5 mg

For Research Use Only

