

## [D - Tyr6, $\beta$ - Ala11, Phe13, Nle14] - Bombesin (6 - 14)

[D - Tyr6,  $\beta$  - Ala11, Phe13, Nle14] - Bombesin (6-14) is a 9 amino acid active fragment of the neuropeptide Bombesin, and is a potent agonist of the orphan bombesin receptor subtype 3 (hBRS-3). Also an agonist of the G-protein-coupled receptors BBR1, BBR2, and BBR3. Bombesin biological effects include hunger control, stimulation of gastrin release from G-cells, hormone release, stimulation of pancreatic enzyme secretion, inhibition of gastric emptying, and modulation of gastric acid secretion.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1158</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 mg per vial
<b>Reconstitution Buffer:</b>	1 mL

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	1298.5 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	$\geq 95\%$
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	$\leq -20^{\circ}\text{C}$ Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	1 mM
<b>Reconstitution Volume:</b>	0.770 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)