

## Glucagon-like Peptide 2 (GLP-2)

Glucagon-like peptide 2 (GLP-2) is a 33 amino acid GI peptide produced upon nutrient consumption via post-translational cleavage of proglucagon in intestinal endocrine cells. It has potent intestinotrophic endocrine/paracrine actions, including intestinal mucosal growth, enhanced brush-border enzyme activity, and delayed gastric transit. May be a potential therapeutic target for subjects with intestinal disease.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1079</b>
<b>Quantity Supplied:</b>	1 vial(s), 0.5 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>	3766.2 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 95% by SDS-PAGE and HPLC
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	≤ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	500 μM
<b>Reconstitution Volume:</b>	0.2655 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)