

β-CGRP, Human

β-calcitonin gene related peptide (β-CGRP) is a potent peptide vasodilator and functions in pain transmission. It is produced in peripheral and central neurons, and when synthesized in the spinal cord may contribute to nervous tissue regeneration after injury.

Materials Provided

Catalog Number:	92-1052
Quantity Supplied:	1 vial(s), 0.5 mg per vial

Description

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

Product Information

Molecular Weight:	3789.4 g/mol
Source:	Synthetic
Purity:	≥ 95% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 μM
Reconstitution Volume:	1.32 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com