

## Bradykinin

Bradykinin is a peptide that acts as a mediator of pain and inflammation, with pro-inflammatory activity. It is an agonist of the bradykinin receptors, with selectivity for B2 receptor over B1. Important roles in fluid regulation, electrolyte balance, smooth muscle contraction, vasodilation and capillary permeability.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1053</b>
<b>Quantity Supplied:</b>	1 vial(s), 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>	1060.2 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 95% by 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>	10 mM
<b>Reconstitution Volume:</b>	0.472 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)