

## Lys-{Des-Arg9}-Bradykinin

Lys-[Des-Arg9]-Bradykinin is a 9 amino acid inflammatory mediating peptide that causes endothelium-dependent vasodilation and hypotensive effects. ACE inhibitors act by increasing bradykinin levels. Potent and selective bradykinin B1 receptor agonist ( $K_i=0.12$  and  $>30000\text{nM}$  for B1 and B2 receptors, respectively). Lys-[Des-Arg9]-bradykinin is 16-fold more potent for B1 than [Des-Arg9]-bradykinin.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1085</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>	1032.2 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	$\geq 95\%$ by HPLC
<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>	0.96 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)