

## WKYMVm-NH2

WKYMVm-NH2 is a peptide found in immune cells that is an agonist of formyl peptide receptors (FPR), with EC50 of 75pM, 75pM, and 3nM for receptors 1, 2, and 3, respectively. It induces calcium mobilization, superoxide production, and chemotactic migration in monocytes and neutrophils. Assists in neutrophil survival and stimulates phospholipase C (PLC) mediated InoP formation.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1075</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 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

<b>Molecular Weight:</b>	856.8 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>	500 µM
<b>Reconstitution Volume:</b>	2.334 mL
<b>Reconstitution Solvent:</b>	H2O

***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)