

## Recombinant Human IL-1 $\alpha$

Interleukin 1, alpha (IL-1 $\alpha$ ) is a pleiotropic cytokine involved in inflammation, immunity, hematopoiesis, and is a pro-apoptotic protein. Stimulates thymocyte proliferation by inducing IL-2 release, B-cell maturation and proliferation, and fibroblast growth factor activity. Has pro-inflammatory properties, also promoting fever and sepsis, and has been implicated in multiple diseases related to these symptoms. Interacts with TNF, another pro-inflammatory cytokine.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1153</b>
<b>Quantity Supplied:</b>	1 vial(s), 2 $\mu$ g 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>	18 kg/mol
<b>Source:</b>	E. coli
<b>Purity:</b>	$\geq$ 98%
<b>Endotoxin Level:</b>	<0.001 ng/mL
<b>Formulation:</b>	
<b>Storage:</b>	$\leq$ -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>	2 $\mu$ M
<b>Reconstitution Volume:</b>	0.055 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)