

## Recombinant Human IL-1 $\beta$

Interleukin-1 beta (IL-1 $\beta$ ) is a secreted proinflammatory cytokine with a broad range of activities including the stimulation of thymocyte proliferation by inducing IL-2 release, B-cell maturation and proliferation, mitogenic FGF-like activity, and the release of prostaglandin and collagenase from synovial cells.

### Materials Provided

|                               |                                |
|-------------------------------|--------------------------------|
| <b>Catalog Number:</b>        | <b>92-1296</b>                 |
| <b>Quantity Supplied:</b>     | 1 vial(s), 10 $\mu$ 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> | 17300 g/mol                                     |
| <b>Source:</b>           | Recombinant                                     |
| <b>Purity:</b>           | $\geq$ 98%                                      |
| <b>Endotoxin Level:</b>  | < 1EU/ $\mu$ g                                  |
| <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>    | 100 $\mu$ g/mL        |
| <b>Reconstitution Volume:</b>  | 0.100 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)