

| Catalog #            | Aliquot Size  |
|----------------------|---------------|
| <b>I860-40N-100</b>  | <b>100 µg</b> |
| <b>I860-40N-250</b>  | <b>250 µg</b> |
| <b>I860-40N-1000</b> | <b>1 mg</b>   |

## Recombinant Human IL1RN

Recombinant human protein expressed in E. coli cells

**Catalog # I860-40N**

Lot # G3211-66

### Product Description

Recombinant Human IL1RN was expressed in E. coli cells. The protein accession number is [P18510](#)

### Gene Aliases

IRAP, ICIL-1RA, IL1 inhibitor, IL1F3, IL-1ra, IL-1RN  
Recombinant Human Interleukin-1 Receptor Antagonist (IL1RN)  
Description

### Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

### Formulation

Lyophilized from a 0.2 µm filtered tris solution with sodium chloride (pH 7.5)

### Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

### Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

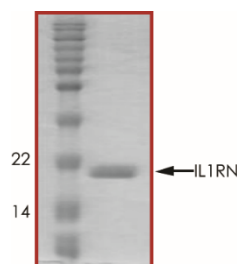
### Scientific Background

IL1RN is found in the conditioned medium of a variety of cells and is produced by macrophages, monocytes, and neutrophils. The soluble form of IL1RN has been shown to be produced by hepatocytes and to be regulated by pro-inflammatory cytokines. It may help in the treatment of sepsis, cachexia, rheumatoid arthritis, chronic myelogenous leukemia, asthma, psoriasis, and inflammatory bowel disease.

### Specific Activity

The ED50 was determined by the ability to inhibit the IL-1 alpha mediated cell proliferation in a murine helper T cell line, for this effect, in the presence of 50 pg/mL of IL-1 alpha, is typically 20-60 ng/mL

### Purity



Recombinant Human IL1RN resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.0 kDa**

## Recombinant Human IL1RN

Recombinant human protein expressed in E. coli cells

|                     |   |
|---------------------|---|
| Catalog Number      | I860-40N  |
| Specific Lot Number | G3211-66  |
| Purity              | 95%   |
| Stability           | 6 mos. at -20°C from date of shipment   |
| Storage & Shipping  | The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**