

## IL1RN Polyclonal Antibody

(Catalog # A52233)

### Background

Inhibits the activity of interleukin-1 by binding to receptor IL1R1 and preventing its association with the coreceptor IL1RAP for signaling. Has no interleukin-1 like activity. Binds functional interleukin-1 receptor IL1R1 with greater affinity than decoy receptor IL1R2; however, the physiological relevance of the latter association is unsure.

### Description

IL1RN Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P18510

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Interleukin-1 receptor antagonist protein (26-177AA)

### Storage

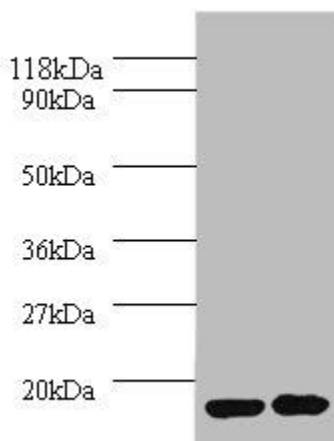
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

ICIL-1RA, IL1 inhibitor, INN=Anakinra IL1RN, IL1F3, IL1RA

### Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



### Western blot

All lanes: IL1RN Polyclonal Antibody at 2ug/ml

Lane 1: EC109 whole cell lysate

Lane 2: 293T whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 21, 18, 20, 17 kDa

Observed band size: 19 kDa