# IL-19 (mouse), (recombinant)

Interleukin-19 (IL-19) is produced by B cells and monocytes. IL-19 binds the interleukin 20 receptor complex to activate STAT3 signaling. IL-19 induces IL-6 and tumor necrosis factor alpha (TNF- $\alpha$ ) expression in monocytes, and promotes T helper 2 cell-mediated immune responses. IL-19 production is up-regulated in resting monocytes following granulocyte-macrophage colony-stimulating factor (GM-CSF) or lipopolysaccharide (LPS) stimulation.

## **Ordering Information**

Order Online »

ENZ-PRT246-0010	10µg
ENZ-PRT246-0100	100µg

Manuals, SDS & CofA

**View Online »** 

### **Handling & Storage**

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For

long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA

solution.

Long Term Storage -80°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Interleukin-19

Endotoxin Content ≤1 EU/μg protein (LAL test)

**Formulation** Lyophilized from a sterile (0.2 micron) filtered solution

containing 10 mM sodium phosphate and 50 mM sodium

chloride, pH 7.5.

**MW** ~17.7 kDa

Purity ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute in sterile water at 0.1 mg/mL.

**Source** Produced in *E. coli*. Contains 153 amino acids.

UniProt ID Q8CJ70

Last modified: May 29, 2024

