# IL-21 (mouse), (recombinant)

Interleukin 21 (IL-21) is normally expressed by activated CD4+ T cells. The IL-21 receptor activates the JAK/STAT signaling pathway and is expressed on T, B, and natural killer (NK) cells. IL-21 cooperates with IL-4 for the production of multiple antibody classes in B cells. IL-21 has pleiotropic effects on the proliferation, differentiation, and effector functions of B, T, NK, and dendritic cells.

## **Ordering Information**

**Order Online** »

ENZ-PRT249-0010	10µg
ENZ-PRT249-0100	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-21

Biological Activity Mouse thymocyte proliferation with costimulation; ED50 ≤

10 ng/mL (≥ 1.0 x  $10^5$  units/mg).

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

Formulation Lyophilized from a sterile (0.2 micron) filtered solution

containing 20 mM sodium bicarbonate, pH 8.5.

**MW** ~15.1 kDa

Purity ≥95% (SDS-PAGE)

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

**Source** Produced in *E. coli*. Contains 130 amino acids.

UniProt ID Q9ES17

Last modified: May 29, 2024



eu@enzolifesciences.com