IL-15 (mouse), (recombinant)

IL-15 is an immunomodulating cytokine that stimulates the proliferation of T lymphocytes and shares many biological properties with IL-2. IL-15 exerts its biological activites primarily on T cells. It is also essential in the development, survival and activation of NK cells. Increased expression of IL-15 has been implicated with rheumatoid arthritis, inflammatory bowel disease and diseases affiliated with retroviruses.

Ordering Information

Order Online »

ENZ-PRT140-0010

10µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-15

Formulation Lyophilized from 10mM sodium phosphate, pH 8.0.

MW ~13.3 kDa

Purity ≥98% (SDS-PAGE gel & HPLC)

Source Produced in *E.coli*.

UniProt ID P48346

Last modified: May 29, 2024

