IL-5 (mouse), (recombinant)

Interleukin 5 (IL-5) is a hematopoietic growth factor that is expressed in type 2 T helper cells, mast cells, and eosinophils. IL-5 acts through the IL-5 receptor, stimulates B cell growth, and mediates eosinophil activation.

Ordering Information

Order Online »

ENZ-PRT240-0025	25µg
ENZ-PRT240-0100	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-5, B-cell differentiation factor I, Eosinophil

differentiaion factor, TFR

Biological Activity TF-1 cell proliferation; ED50 \leq 250 ng/mL (\geq 4.0 x 10⁶

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 9.

MW Dimer, ~13.1/26.3 kDa

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute in sterile water at 0.1 mg/mL.

Source Produced in *E. coli*. Contains 113/226 amino acids.

UniProt ID P04401

Last modified: May 29, 2024

