IL-17A (mouse), (recombinant)

Interleukin 17A (IL-17A), also known as CTLA-8, is a member of the IL-17 family of proteins. IL-17A is a pro-inflammatory cytokine that is secreted by activated CD4+ and CD8+ T lymphocytes. IL-17A acts through its receptor, IL-17R, to promote increased cytokine and chemokine secretion. Cytokines and chemokines then mediate the immunoregulatory function of IL-17A by promoting the proliferation, maturation, and chemoattraction of neutrophils to inflammatory sites. Elevated levels of IL-17A are associated with rheumatoid arthritis, airway inflammation, allograft rejection, inflammatory bowel disease, psoriasis, cancer, and multiple sclerosis.

Ordering Information

Order Online »

ENZ-PRT244-0025	25µg
ENZ-PRT244-0100	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-17A, IL-17, CTLA-8, Cytotoxic T-lymphocyte-

associated antigen 8

Biological Activity Production of IL-6 by mouse 3T3 cells; ED50≤5 ng/mL

(≥2.0×10⁵ units/mg).

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

Formulation Lyophilized from a sterile (0.2 micron) filtered solution

containing 0.1% Trifluoroacetic Acid (TFA).

MW Dimer, ~15.1/30.2 kDa

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute in sterile water at 0.1 mg/mL.

Source Produced in *E. coli*. Contains 134/268 amino acids.

UniProt ID Q62386

Last modified: May 29, 2024

