

Human IL-17A (recombinant)



Alias: none **Catalog #:** 6587

Size: 5 ug **Research Use Only**

Molecular Weight: 15.1 kDa

Source: Yeast. Recombinant Human IL-17A was produced in yeast and, therefore, does not have endotoxin. It is naturally folded and post-translationally modified.

Formulation: Lyophilized without carrier protein.

Purity: >95% as visualized by SDS-PAGE analysis.

Purification: Ion-exchange chromatography.

Bioactivity: In testing

Entrez Gene ID: 3605

Number of Amino Acids: 132

Amino Acid Sequence: GITIPRNPGC PNSDKNFPR TVMVNLNIHN RNTNTNPKRS SDYYNRSTSP WNLHRNEDPE
RYPSSIWEAK CRHLGCINAD GNVDYHMNSV PIQQEILVLR REPPHCPNSF RLEKILVSVG
CTCVTPIVHH VA (132)

Country of Origin: USA

Reconstitution: Reconstitute with sterile phosphate-buffered saline containing at least 0.1% carrier protein.

Stability and Storage: Stable for up to twelve months from date of receipt at -20°C. Stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C. Avoid repeated freeze/thaw cycles.

Applications: The human IL-17A protein can be used in cell culture, as a IL-17A ELISA Standard, and as a Western Blot Control.

Background: IL-17A is a member of the IL-17 family, which is comprised of 6 members [IL-17A, IL-17B, IL-17C, IL-17D, IL-17E (also called IL-25), and IL-17F]. IL-17 family members are involved in numerous immune regulatory functions, including inducing and mediating proinflammatory responses and allergic responses. IL-17 induces the production of many other cytokines (IL-6, G-CSF, GM-CSF, IL-1beta, TGF-beta, and TNF-alpha), chemokines, including IL-8 (CXCL8), GRO-alpha (CXCL1) and MCP-1 (CCL2) and prostaglandins from many cell types (fibroblasts, endothelial cells, epithelial cells, keratinocytes and macrophages).

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Human IL-17A updated on: 8/21/2014