

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.

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 (≥2.0×10 ⁵ 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 <i>E. coli</i> . Contains 134/268 amino acids.
UniProt ID	Q62386

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)