

IL-17A (human), (recombinant)

IL-17A is produced as a disulfide-linked homodimer that can also form heterodimers with IL-17F. IL-17 acts via ubiquitously expressed IL-17 receptors, inducing expression of pro-inflammatory factors such as IL-1, IL-6, TNF, CXCL8/IL-8, and matrix metalloproteinases. Elevated levels of IL-17A and increased presence of Th17 cells has been associated with airway inflammation, rheumatoid arthritis, intraperitoneal abscesses, inflammatory bowel disease, allograft rejection, psoriasis, multiple sclerosis, and cancer.

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-17, CTLA8, Cytotoxic T-lymphocyte-associated serine esterase 8

Formulation Lyophilized with no additives.

MW ~31.3 kDa

Purity ≥98% (SDS & HPLC)

Purity Detail Sterile filtered through a 0.2 micron filter.

Source Produced in *E. coli*.

UniProt ID Q16552

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)