

# 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.

## Ordering Information

[Order Online »](#)

ENZ-PRT188-0025	25µg
-----------------	------

## 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



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)