

# IL-13 (human), (recombinant)

IL-13 (interleukin-13) is primarily produced by activated T cells and has many of the same biological properties as IL-4. IL-13 suppresses the expression of pro-inflammatory cytokines such as IL-1, IL-6, TNF- $\alpha$  and IL-8 by monocytes/macrophages, induces CD23 expression on B cells, promotes B cell proliferation, stimulates secretion of IgE and IgG4, and induces IFN- $\gamma$  production by NK cells. IL-13 has been shown to inhibit the replication of HIV both in vitro and in vivo. IL-13 induces many phenotypes of allergic lung diseases, including airway hyper responsiveness, goblet cell metaplasia and mucus hypersecretion that contribute to airway obstruction.

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-PRT183-0010	10 $\mu$ g
-----------------	------------

## 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-13, NC30

**Biological Activity** Determined by its ability to stimulate the proliferation of human TF-1 cells. The expected ED50 is  $\leq 1.0$  ng/ml, corresponding to a specific activity of  $\geq 1 \times 10^6$  units/mg.

**Endotoxin Content** Endotoxin level is  $< 0.1$  ng/ $\mu$ g of protein ( $< 1$  EU/ $\mu$ g)

**Formulation** Lyophilized from 20mM sodium phosphate, pH 7.0.

**MW** ~12.6 kDa

**Purity**  $\geq 98\%$  (SDS & HPLC)

**Purity Detail** Sterile filtered through a 0.2 micron filter.

**Source** Produced in *E. coli*.

**UniProt ID** P35225



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)