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-α 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-γ 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

10µg

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 x 10⁶ units/mg.

Endotoxin Content Endotoxin level is < 0.1 ng/μg of protein (< 1 EU/μg)

Formulation Lyophilized from 20mM sodium phosphate, pH 7.0.

MW ~12.6 kDa

Purity ≥98% (SDS & HPLC)

Purity Detail Sterile filtered through a 0.2 micron filter.

Source Produced in E. coli.

UniProt ID P35225

Last modified: May 29, 2024

