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

Last modified: May 29, 2024



uk@enzolifesciences.com