

Catalog #	Aliquot Size
1896-40N-20	20 µg
1896-40N-100	100 µg
1896-40N-1000	1 mg

Recombinant Human IL36A

Recombinant human protein expressed in E. coli cells

Catalog # 1896-40N

Lot # G3211-72

Product Description

Recombinant Human IL36A was expressed in E. coli cells. The protein accession number is [Q9UHA7](#)

Gene Aliases

IL-1 Epsilon, FIL1 epsilon, Interleukin-1 family member 6, IL-1F6, IL-36 gamma, IL-1F, IL-1H1, Recombinant Human Interleukin-36 alpha (IL36A)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

Formulation

Lyophilized from a 0.2 µm filtered solution in PBS, NaCl and Tween® 20 with BSA as a carrier protein

Reconstitution Protocol

Reconstitute at 10 µg/mL in sterile PBS containing at least 0.1% human or bovine serum albumin

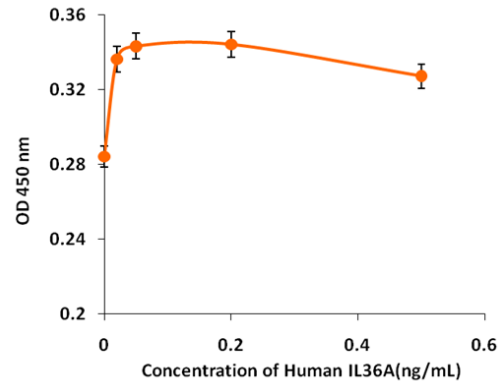
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

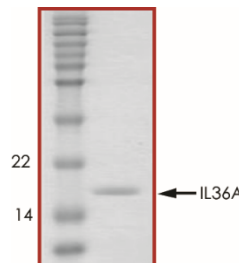
Scientific Background

IL-1 epsilon is an actively secreted member of the IL-1 family of cytokines. It is mainly expressed in tissues with epithelial cells and in psoriatic plaques, as well as by monocytes, B cells and T cells. IL-1 epsilon has been shown to signal through the IL-Rrp2 receptor which is primarily expressed on certain dendritic cells, and to act as an agonist of NF-kappaB which then stimulates the inflammatory response in bronchial epithelial cells.

Specific Activity



Purity



Recombinant Human IL36A resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.4 kDa**

Recombinant Human IL36A

Recombinant human protein expressed in E. coli cells

Catalog Number	1896-40N
Specific Lot Number	G3211-72
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.