

Recombinant Human IL36B

Recombinant human protein expressed in E. coli cells

Catalog # 1896-40BN

Lot # G3211-73

Product Description

Recombinant Human IL36B was expressed in E. coli cells. The protein accession number is [Q9NZH7](#)

Gene Aliases

IL1 eta, Interleukin-1 eta, Interleukin-1 family member 8, Interleukin-1 homolog 2

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 (with BSA as a carrier protein)

Reconstitution Protocol

Reconstitute in sterile PBS at 100µg/mL (containing at least 0.1% human or bovine serum albumin)

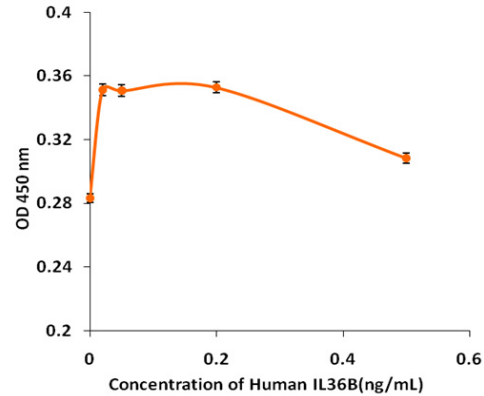
Storage and Stability

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

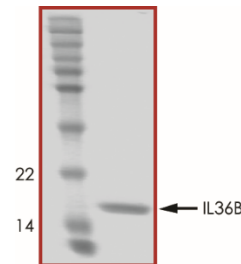
IL36B, also known as Interleukin 36 beta and IL1F8, is a cytokine belonging to the IL-1 family. It is actively secreted and is found to form a cytokine gene cluster on chromosome 2 with eight other IL1 cytokines. IL36B signals through the IL1Rrp2/IL-36R receptor on dendritic cells which then activates NFκB and MAPK signalling pathways to create a pro-inflammatory response. It stimulates the production of mature adipocytes, and IL6 and IL8 synovial fibroblasts. It also induces expression of some antimicrobial peptides and matrix metalloproteases. It is highly expressed in psoriatic plaques, as part of the IL36 signalling system is found in many epithelial barriers.

Specific Activity



IL-36 beta can bind to human IL-1 Rrp2 Fc Chimera with a linear range of 0.15-5 µg/mL as determined by its binding ability in a functional ELISA

Purity



Recombinant Human IL36B resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.5 kDa**

Recombinant Human IL36B

Recombinant human protein expressed in E. coli cells

Catalog Number	1896-40BN
Specific Lot Number	G3211-73
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.