

## Recombinant Human IL17F

Recombinant human protein expressed in E. coli cells

**Catalog # I877-40FN**

Lot # G3211-61

### Product Description

Recombinant Human IL17F was expressed in E. coli cells. The protein accession number is [Q96PD4](#)

### Gene Aliases

Cytokine ML-1

### 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 Acetonitrile and TFA (with BSA as a carrier protein)

### Reconstitution Protocol

Reconstitute in sterile 4mM HCl at 25µg/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.

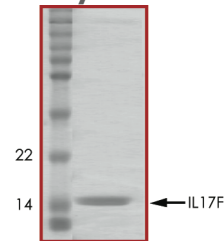
### Scientific Background

Interleukin 17F is a disulfide-linked homodimeric protein belonging to the IL-17 family of proinflammatory cytokines. Members belonging to this family share conserved features of cysteine-knot structures but differ in their N-terminal regions. The T-helper cell produced IL-17 family is inducible by IL15 and IL23, and IL17F induces T-cell and PBMC activation and proliferation. It also performs cartilage matrix turnover and angiogenesis prevention. Out of the IL-17 family, IL17F is the most homologous to IL17 and dimerizes in a way similar to that of nerve growth factors (NGF) and other neurotrophins.

### Specific Activity

The ED50 is 8.0-40.0ng/mL as determined by its ability to induce IL-6 secretion by NIH-3T3 mouse embryonic fibroblast cells.

### Purity



Recombinant Human IL17D resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **15.2 kDa**

## Recombinant Human IL17F

Recombinant human protein expressed in E. coli cells

Catalog Number	I877-40FN
Specific Lot Number	G3211-61
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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