

## Recombinant Human WISP3

Recombinant human protein expressed in E. coli cells

**Catalog # W810-40N**

Lot # G3223-28

### Product Description

Recombinant Human WISP3 was expressed in E. coli cells. The protein accession number is [O95389](#)

### Endotoxin Level

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

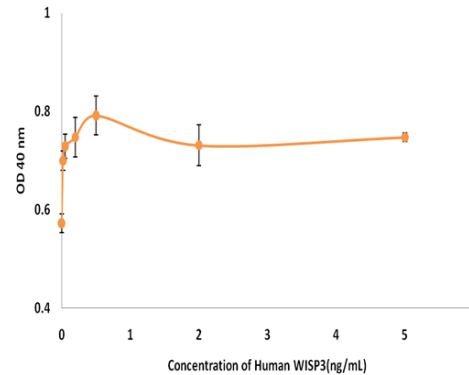
### Storage and Stability

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

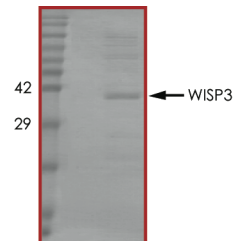
WNT1-inducible-signaling pathway protein 3 (WISP3) is a member of the CCN family of cysteine-rich, glycosylated signalling proteins that mediate several developmental processes, and a member of the WNT1 inducible signalling pathway (WISP) subfamily. Members of this family are characterized by four conserved cysteine-rich domains: insulin-like growth factor-binding domain (IGFBP), von Willebrand factor type C (VWFC) domain, thrombospondin type-1 repeat (TSP type-1) domain, and C-terminal cystine knot-like (CTCK) domain. WISP3 is expressed in mature and fetal kidneys, testis, and in lower levels in the ovary, placenta, prostate, and small intestine. It is necessary for the growth of postnatal skeleton and cartilage homeostasis.

### Specific Activity



The ED50 was 0.2-0.3 µg/mL as determined by cell proliferation assay

### Purity



Recombinant Human WISP3 resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **37.1 kDa**

## Recombinant Human WISP3

Recombinant human protein expressed in E. coli cells

Catalog Number

**W810-40N**

Specific Lot Number

**G3223-28**

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.**