



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Human Oncostatin M

Catalog #: 90227-B (10 µg)
Lot #: 150729

Description: Recombinant Oncostatin-M is a disulfide-linked homodimeric protein consisting of 228 amino acid residues, and migrates as an approximately 26 kDa protein under non-reducing and reducing conditions in SDS-PAGE. Optimized DNA sequence encoding Human Oncostatin-M mature chain was expressed in *E. coli*.

Source: Optimized DNA sequence encoding human Oncostatin M mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 µm filtered concentrated (1 mg/ml) PBS solution, pH 7.0.

Reconstitution: Reconstitute at 0.1-1.0 mg/ml in distilled water. This solution can then be diluted into other buffers. To maximize product collection from vial surface, vortex briefly and then spin down to recollect the liquid.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at +4°C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: ≥95% by SDS-PAGE and HPLC

Endotoxin Level: <0.1 ng/µg (1 EU/µg), using the LAL gel clot method.

Biological Activity: The ED₅₀ was determined by the dose-dependent stimulation of the proliferation of human TF-1 cells is ≤3 ng/ml, corresponding to a specific activity of ≥2x10⁵ units/mg.

Protein Sequence:

```
AAIGSCSKEYRVLLGQLQKQTDLMQD  
TSRLDPYIRIQGLDVPKLRHCRERP  
GAFPSEETLRGLGRRGFLQTLNATLG  
CVLHRLADLEQRLPKAQDLERSGLNIE  
DLEKLMARPNIPLGRNNIYCMAQLLD  
NSDTAEPTKAGRGASQPPTPTPASDA  
FQRKLEGCRFLHGYHRFMHSVGRVFS  
KWGESPNRSRRHSPHQALRKGVRRT  
RPSRKGRMLMTRGQLPR
```

References:

1. Am. J. Respir. Cell Mol. Biol., May 2009, 40: 620 – 630
2. J. Virol., Apr 2009, 83: 3298 - 3311.
3. Clin. Cancer Res., Mar 2009, 15: 1519 - 1526.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com