



6405 Mira Mesa Blvd., Suite 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

## IL-11, Human

Catalog #: 90176-B  
Lot#: 220621

**Description:** Recombinant human interleukin-11 (IL-11), also known as oprelvekin, is a disulfide-linked monomeric protein consisting of 180 amino acids and migrates as an approximately 19 kDa protein under reducing and non-reducing conditions.

**Source:** Optimized DNA sequence encoding human IL-11 mature chain was expressed in *E. coli*.

**Formulation:** Lyophilized from 0.2 µm filtered phosphate buffered solution in 2.5% glycine, pH 4.5.

**Reconstitution:** A quick spin of the vial followed by reconstitution in distilled water to concentration not less than 0.1 mg/ml. This can then be diluted into other buffers.

**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 2 - 8°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:** >97%, as determined by SDS-PAGE and HPLC.

**Endotoxin Level:** Endotoxin level was found to be < 0.1 ng/µg (1EU/µg), using the LAL gel clot method.

**Biological Activity:** The ED50, determined by the dose-dependent stimulation of the proliferation of murine T11 cells is ≤1.0 ng/ml, corresponding to a specific activity of ≥1 x 10<sup>6</sup> units/mg.

**Protein Sequence:**  
PGPPPGPPRVSPDPRAELDSTVLLTR  
SLLADTRQLAAQLRDKFPADGDHNL  
SLPTLAMSAGALGALQLPGVLTRLRAD  
LLSYLRHVQWLRRRAGGSSSLKTLEPEL  
GTLQARLDRLRLQLLMSRLALPQPP  
PDPPAPPLAPPSSAWGGIRAHAAILGG  
LHLTLDWAVRGLLLLKTRL

### References:

1. Dimitriadis, E., *et al.* (2010). *Placenta* **31 Suppl**:S99-104.
2. Nagineni, C.N., *et al.* (2010). *Biochem. Biophys. Res. Commun.* **391(1)**:287-92.