



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-A

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 10⁶ units/mg.

Protein Sequence:

PGPPPGPPRVSPDPRAELDSTVLLTR
SLLADTRQLAAQLRDKFPADGDHNLD
SLPTLAMSAGALGALQLPGVLRRLRAD
LLSYLRHVQWLRRLRAGGSSSLKTLEPEL
GTLQARLDRLRLRLQLLMSRLALPQPP
PDPPAPPLAPPSSAWGGGIRAAHAILGG
LHLTLDWAVRGLLLLLKTRL

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.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com