



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

## Rat Interleukin 10 Recombinant

Catalog #: 90175-A  
Lot#: 230817

**Description:** Recombinant Rat Interleukin-10 is a disulfide-linked monomer protein consisting of 161 amino acid residues, and migrates as an approximately 19 kDa protein under non-reducing and reducing conditions.

**Source:** Optimized DNA sequence encoding rat IL-10 mature chain was expressed in *E. coli*.

**Formulation:** Lyophilized from 0.2 µm filtered PBS.

**Reconstitution:** A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution 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:** >95%, 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 inhibition of antigen-specific T cell proliferation, is less than 12 ng/ml.

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](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)