



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

Mouse Interleukin-1 beta Recombinant

Catalog #: 90172-B
Lot#: 230131

Description: Recombinant murine interleukin-1 β (IL-1 β), also known as catabolin, is a disulfide-linked monomeric protein consisting of two 153 amino acid residue subunits and migrates as an approximately 17 kDa protein under reducing and non-reducing conditions.

Source: Optimized DNA sequence encoding murine IL-1 β mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 μ m filtered solution in PBS, pH 7.0.

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: >98%, as determined by SDS-PAGE and HPLC.

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

Biological Activity: The ED₅₀, determined by the dose-dependent stimulation of thymidine uptake by murine D10S cells, is \leq 2.0 pg/ml, corresponding to a specific activity of \geq 5 x 10⁸ units/mg.

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