



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

IL-2RB Recombinant (CD122)

Catalog #: 79655
Lot#: 181023

Description: Recombinant IL2RB is a monomer protein consisting of 451 amino acid residue subunits, due to glycosylation migrates as an approximately 60kDa protein on SDS-PAGE.

Source: Optimized DNA sequence encoding extracellular domain of human Interleukin-2 receptor beta subunit including an Fc tag was expressed in HEK293 cells.

Formulation: Lyophilized from a 0.2 µm filtered PBS solution.

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 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: >90%, 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: Measured by its ability to bind biotinylated recombinant rat IL2 in a functional ELISA.

References:

1. Zheng, C. *J. Immunol.* Jun 2008; **180**: 7221-7229.
2. Jin, H. *Int. Immunol.*, Apr 2006; **18**: 565 - 572.