

Description: Recombinant human Interleukin-6, also known as IL-6, is a disulfide-linked homodimeric protein consisting of 183 amino acids and migrates as an approximately 20 kDa protein under reducing conditions.

Source: Optimized DNA sequence encoding human IL-6 mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 µm filtered PBS solution with 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: >97%, as determined by SDS-PAGE and HPLC.

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

Biological Activity: The ED50, determined by the dose-dependent stimulation of the proliferation of murine 7TD1 cells was found to be less than 0.1 ng/ml, corresponding to specific activity of 1×10^7 IU/mg.

References:

1. *Glycobiology*, Oct 2009; 19: 1082 - 1093.
2. *Lab Med*, Oct 2009; 40: 600 - 603.