



GIBCO BRL

LIFE  TECHNOLOGIES™

Recombinant Human Interleukin-2 (IL-2)

Cat. No.: 13238-050

Size: 5 µg

Lot No.:

Store at -20°C.

Description:

Recombinant Human Interleukin-2, from *E. coli*, is a 15-kDa protein containing 133 amino-acid residues (1). It is a potent lymphoid cell growth factor exerting its biological activity on T lymphocytes, large granular lymphocytes, monocytes, B lymphocytes and oligodendritic cells (2,3). For most *in vitro* applications, IL-2 exerts its biological activity in the concentration range of 0.1 to 1 ng/ml. Recombinant Human IL-2 is provided as a lyophilized protein obtained from a membrane-filtered (0.22 µM) solution of distilled water with no additives.

Directions for Use:

1. Reconstitute in 100 mM acetic acid to a concentration of 100 ng/µl. **Note:** **Do not store in glass;** store in polypropylene vial. Store at -20°C reconstituted for up to six months. Avoid repeated freeze-thaw cycles.
2. Reconstituted solution can be further diluted in distilled water or aqueous buffers to working concentration. **Do not store dilute solutions.**

Note: To avoid adsorptive loss, dilute solutions should be supplemented with Bovine Serum Albumin (BSA) at a concentration of not less than 0.1%.

Doc. Rev.: 080197

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE™ (800) 828-6666.

Quality Control:

- Purity: $\geq 97\%$ (by SDS-PAGE)
- Endotoxin: ≤ 0.1 ng/ μ g of IL-2
- Specific Activity: units/mg
- Biological Activity: $ED_{50} \leq 0.1$ ng/ml (determined by dose-dependent stimulation of murine CTLL-2 cells)

References:

1. Morgan, D.A. *et al.* (1976) *Science* 193, 1007.
2. Ruscetti, F.W. *et al.* (1977) *J. Immunol.* 119, 131.
3. Lotze, M.T. (1992) *Interleukin-2 In: Human Cytokines*, Eds. B.B. Aggarwal, and J.U. Gutterman, Blackwell Scientific Press, Boston, MA, p.81.