

IL-4

**Interleukin 4, B-cell stimulatory factor
human, recombinant, *E. coli***

Cat. No.	Amount
PR-463	10 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Lyophilized.

Solubility

It is recommended to reconstitute the lyophilized IL-4 in bidest H₂O not less than 100 µg/ml, which can then be further diluted to other aqueous solutions. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

Activity

ED₅₀: < 0.2 ng/ml, corresponding to a specific activity of 5 × 10⁵ IU/mg, determined by the dose-dependent stimulation of TF-1 cells.

Purity

≥ 98% by SDS-PAGE, RP-HPLC, and FPLC.

Endotoxin

Less than 0.1 ng/µg (IEU/µg) of IL-4.

Description

The anti-inflammatory cytokine Interleukin-4 is involved in the regulation of inflammatory responses by promoting the differentiation of naïve T-cells to T helper type 2 (Th2) cells.

A key anti-inflammatory action of IL-4 results from its ability to inhibit the release of pro-inflammatory cytokines by innate immune cells and to upregulate the synthesis of IL-1 receptor antagonist.

IL-4 has been demonstrated to be a potent cofactor for B and T lymphocyte proliferation and differentiation.

Recombinant human IL-4 produced in *E. coli* is a single, non-glycosylated form of human IL-4 polypeptide chain containing 130 amino acids and having a molecular mass of 15 kDa.

IL-4 is purified by proprietary chromatographic techniques.

Selected References:

- Wurster *et al.* (2002) Interleukin-4-mediated protection of primary B cells from apoptosis through Stat6-dependent up-regulation of BclxL. *J. Biol. Chem.* **277**:27169.
- Nolan *et al.* (2004) Role of interleukin-4 in regulation of age-related inflammatory changes in the hippocampus. *J. Biol. Chem.* **280**:9354.
- Petersen *et al.* (2002) Optimisation of the CT h4S bioassay for detection of human interleukin-4 secreted by mononuclear cells stimulated by phytohaemagglutinin or by human leukocyte antigen mismatched mixed lymphocyte culture. *Immunol. Lett.* **84**:29.
- Whitehead *et al.* (2002) Phase II trial of recombinant human interleukin-4 in patients with advanced renal cell carcinoma: a southwest oncology group study. *J. Immunother.* **25**:352.
- Ekerfelt C. and Ernerudh J. (2002) Detection of spontaneous and antigen-induced human interleukin-4 responses in vitro: comparison of ELISPOT, a novel ELISA and real-time RT-PCR. *J. Immunol. Methods.* **260**:55.
- Casolaro *et al.* (2000) Identification and characterization of a critical CP2-binding element in the human interleukin-4 promoter. *J. Biol. Chem.* **275**:36605.
- Davis *et al.* (2000) A phase I and pharmacokinetic study of subcutaneously-administered recombinant human interleukin-4 (rhIL-4) in patients with advanced cancer. *Growth Factors.* **17**:287.
- Okada *et al.* (2000) Characterization and transduction of a retroviral vector encoding human interleukin-4 and herpes simplex virus thymidine kinase for glioma tumor vaccine therapy. *Cancer Gene Ther.* **7**:486.