

Interleukin-2, human (hIL-2)

recombinant (*E. coli*)

Solution, filtered through 0.2 µm pore size membrane

Version 17
Content version: February 2019
Store at -15 to -25°C

Cat. No. 10 799 068 001 10,000 U (5 µg, 50 ml)
Cat. No. 11 011 456 001 10,000 U (5 µg, 1 ml)
Cat. No. 11 147 528 001 50,000 U (25 µg, 5 ml)

1. What this Product Does

Contents

Cat. No.	U	µg	ml	U/ml	µg/ml
10 799 068 001	10,000	5	50	200	0.1
11 011 456 001	10,000	5	1	10,000	5
11 147 528 001	50,000	25	5	10,000	5

Solution in PBS (phosphate buffered saline) and 1 mg/ml BSA (bovine serum albumin), [purity of BSA: >98%, endotoxin (LAL): <1 EU/mg BSA], filtered through 0.2 µm pore size membrane.

Storage and Stability

Stable at -15 to -25°C until the expiration date printed on the label. Store the solution in aliquots at -15 to -25°C.
⊗ Avoid repeated freezing and thawing.

Application

Recombinant IL-2, human allows

- the cultivation of human and murine IL-2 dependent T-cell lines and natural killer cell lines,
- the proliferation of mitogen-activated T-lymphocytes and natural killer cells,
- the establishment of human and murine thymocyte, splenocyte or peripheral blood lymphocyte (PBL) derived T-cell lines, and
- the generation of human and murine lymphokine-activated killer (LAK) cells (1-11).

2. How to Use this Product

2.1 Before you Begin

Working Concentration

Established IL-2-dependant T-cell lines usually require 10-20 U/ml. Add IL-2 to the freezing medium for IL-2 dependant cell lines.

Recommended Method of Dilution

Dilute the concentrated IL-2 solution (200 U/ml or 10,000 U/ml) with PBS or culture medium containing 1 mg/ml BSA or HSA (0.1%) or 1 - 10% serum.

Reagents Required

- Culture medium, e.g. RPMI 1640, containing 10% FCS, 10 mM Hepes*, 2 mM L-glutamine, (1×) non-essential amino acids, 1 mM sodium pyruvate and 50 mM β-mercaptoethanol.
- Cell Proliferation Kit II (XTT)*

2.2 Procedure

Instructions for the Determination of hIL-2 Activity on IL-2 Dependent Cells (IL-2 Proliferation Assay; Exemplary with CTLL-2 cells)

- 1 Prepare hIL-2 test samples as serial dilutions (final concentration e.g., from 0.001 - 100 ng/ml) in culture medium in a 96-well flat-bottomed microplate (tissue culture grade) in a volume of 100 µl.
⊗ As a negative control use culture medium alone.
- 2 Harvest sensitive cells, e.g., CTLL-2 cells (ATCC TIB 214) and wash them three times by centrifugation in culture medium without hIL-2. Resuspend and adjust the cells in culture medium to 4×10^4 cells/ml.
- 3 Add 100 µl of the cell suspension to 100 µl of the prediluted hIL-2 samples into each well, revealing a final cell number of 2×10^4 cells/ml (4×10^3 cells/well).
- 4 Incubate the microplate for 48 h at 37°C and 5% CO₂.
- 5 After the incubation period add 100 µl XTT labeling mixture [for details see working instruction of the Cell proliferation kit II (XTT)] to each well. Incubate the microplate for another 6 h at 37°C and 5% CO₂.
- 6 Measure the spectrophotometrical absorbance of the samples by using a microplate (ELISA) reader. To measure absorbance of the formazan product a wavelength between 450 and 500 nm (e.g., 492 nm) can be used, depending on the optical filters available for the ELISA reader used. The reference wavelength should be more than 650 nm.

3. Results

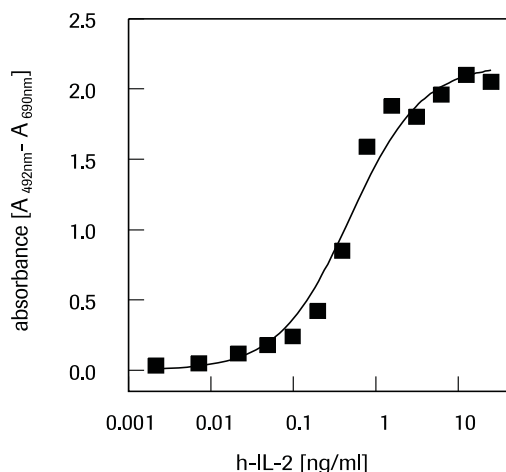


Fig. 1: Stimulation of cell proliferation with mouse CTLL-2 cells in response to recombinant IL-2, human (h-IL-2) using the procedure described.

4. Additional Information on this Product

Product Description

Interleukin-2 (IL-2, also known as T-Cell Growth Factor, TCGF) is a lymphokine which is produced by lectin- or antigen-activated T-cells and plays an important immunoregulatory role. This factor, or lymphokine, was first identified by its ability to promote the long-term *in vitro* proliferation of activated T-cells. It also promotes the generation and proliferation of cytotoxic T-cells, natural killer (NK) cells, and lymphokine-activated killer (LAK) cells (1-11). Recombinant human IL-2 allows the cultivation of human and murine IL-2-dependent T-cell lines and natural killer cell lines, the proliferation of mitogenactivated T-lymphocytes and natural killer cells, the establishment of human and murine thymocyte-, splenocyte-, or peripheral blood lymphocyte (PBL)- derived T-cell lines, and the generation of human and murine lymphokine-activated killer (LAK) cells (1-11).

Preparation

Recombinant interleukin-2, human (hIL-2) is produced in *E. coli* and purified by standard chromatographic techniques.

Primary Structure

The primary structure of recombinant, IL-2, human is identical to that of natural, human IL-2 (one polypeptide chain, 133 amino acids), but recombinant IL-2 has an extra methionine at the amino-terminus (one polypeptide chain, 134 amino acids) and is not glycosylated (12-18). Glycosylation is not essential for biological activity.

Purity

≥95% pure as determined by SDS-PAGE [Endotoxin level: <0.1 EU/μg (LAL-test), <10 EU/ml (LAL-test)].

⊙ 1 EU corresponds to 0.1 ng.

Specific Activity/EC₅₀

>2 × 10⁶ U/mg (<0.5 ng/ml) [(hIL-2, NIBSC, 1st international standard, 86/504), at least the same specific activity (EC₅₀) compared to the indicated standard is guaranteed](19,20). Human, recombinant IL-2 has the same biological activity *in vitro* as compared to human, natural IL-2 (15, 1-4).

E₅₀ Definition/Unit Definition

The amount of hIL-2 that is required to support half-maximal stimulation of cell proliferation (XTT cleavage) with CTLL-2 cells (1 unit equals ≤0.5 ng).

Molecular Weight

15,000 Da

Species Specificity

Recombinant IL-2, human is effective on mouse and human cells.

References

- 1 Talmadge, J. E. (1985) *J. Biol. Resp. Modif.* **4**, 18-34.
- 2 Naruo, K. et al. (1985) *Biochem. Biophys. Res. Commun.* **128**, 257.
- 3 Rosenberg, S. A. et al. (1984) *Science* **223**, 141.
- 4 Roifman, C. M. et al. (1985) *Cell. Immunol.* **95**, 14.
- 5 Gillis, S. (1983) *J. Clin. Immunol.* **3**, 1-13.
- 6 Smith, K. A. (1984) *Ann. Rev. Immunol.* **2**, 319-333.
- 7 Cantrell, D. A. & Smith, K. A. (1984) *Science* **224**, 1312-1316.
- 8 Fletcher, M. & Goldstein, A. L. (1987) *Lymphokine Res.* **6**, 43-57.
- 9 Smith, K. A. (1988) *Science* **240**, 1169-1176.
- 10 Hatakeyama, M. & Taniguchi, T. (1990) In: Peptide Growth Factors and Their Receptors I (Sporn, M. B. & Roberts, A. B., eds.) Springer Verlag Berlin Heidelberg New York, pp. 523-540.
- 11 Winkelhake, J. L. & Gaung, S. S. (1990) *Pharmacol. Rev.* **42**, 1-28.
- 12 Devos, R. et al. (1983) *Nucleic Acids Res.* **11**, 4307.
- 13 Taniguchi, T. et al. (1983) *Nature* **302**, 305-310.

- 14 Maeda, S. et al. (1983) *Biochem. Biophys. Res. Commun.* **115**, 1040-1047.
- 15 Liang, S.-M. et al. (1985) *Biochem. J.* **229**, 429-439.
- 16 Fujita, T. et al. (1983) *Proc. Natl. Acad. Sci. USA*, **80**, 7437-7441.
- 17 Degraeve, W. et al. (1983) *EMBO J.* **2**, 2349-2353.
- 18 Robb, R. J. (1985) *Methods Enzymol.* **116**, 493-525.
- 19 Gearing, A. J. H., Johnstone, A. P. & Thorpe, R. (1985) *J. Immunol. Methods* **83**, 1.
- 20 Gearing, A. J. H. & Thorpe, R. (1988) *J. Immunol. Meth.* **114**, 3-9.

* available from Roche Diagnostics

5. Supplementary Information

Changes to Previous Version

- Editorial changes.

Text Conventions

To make information consistent and understandable, the following text conventions are used in this Instruction Manual:

Text Convention	Use
Numbered instructions labeled 1 , 2 etc.	Steps in a procedure that must be performed in the order listed.

Symbols

Symbols are used in this Instruction Manual to highlight important information:

Symbol	Description
⊙	Information Note: Additional information about the current topic or procedure.

Trademarks

All third party product names and trademarks are the property of their respective owners.

Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit sigma-aldrich.com, and select your home country. Country-specific contact information will be displayed.



Roche Diagnostics GmbH
Sandhofer Strasse 116
68305 Mannheim
Germany