

# Interleukin-1 $\beta$ , human (hIL-1 $\beta$ )

recombinant (*E. coli*)

Solution, filtered through 0,2  $\mu$ m pore size membrane

Cat. No. 11 457 756 001

100,000 U (2  $\mu$ g, 1 ml)

**Version 04**  
Content version: June 2011  
Store at -15 to -25°C

## 1. What this Product Does

### Contents

100,000 U/ml (2  $\mu$ g/ml) 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  $\mu$ m 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

Interleukin-1 (IL-1), secreted by activated monocytes or macrophages and other cell types, is a pleiotropic factor for a variety of sensitive cell.

## 2. How to Use this Product

### 2.1 Before you Begin

#### Working Concentration

hIL-1 $\beta$  exerts its biological activity in the concentration range of 5–500 U/ml (0.1–1 ng/ml).

#### Recommended Method of Dilution

Dilute the concentrated IL-1 $\beta$  solution (100,000 U/ml) with PBS or culture medium containing 1 mg/ml BSA [or HSA (human serum albumin)] or 1–10% serum.

#### Reagents Required

- Culture medium, e.g. RPMI 1640, containing heat inactivated 10% FCS, 10 mM HEPES\*, 2 mM L-glutamine, (1 $\times$ ) non-essential amino acids, 1 mM sodium pyruvate and 50  $\mu$ M  $\beta$ -mercaptoethanol and 50 U/ml interleukin-2\*.
- hIL-1 $\beta$  stock solution (100,000 U/ml; 2  $\mu$ g/ml); [ $^3$ H]-thymidine.

### 2.2 Procedure

#### Instructions to determine the activity of recombinant human IL-1 $\beta$ on sensitive cells (with mouse C3H/HeJ thymocytes as an example)

- 1 Seed mouse C3H/HeJ thymocytes at a concentration of  $5 \times 10^5$  cells/well in 200  $\mu$ l culture medium containing various amounts of hIL-1 $\beta$  [final concentration 0.5–500 U/ml (0.01–10 ng/ml)] into microplates (tissue culture grade, 96 wells) and incubate for 3 days at 37°C and 6.5% CO<sub>2</sub> in a humidified atmosphere.
- 2 Add 1  $\mu$ Ci/well [ $^3$ H]-thymidine, and incubate for another 24 h as described.
- 3 After this incubation period, harvest the contents of each well onto glass fiber filters using a cell harvester, and determine the radioactivity incorporated into the DNA using an  $\alpha$ - $\beta$ -scintillation counter.

## 3. Results

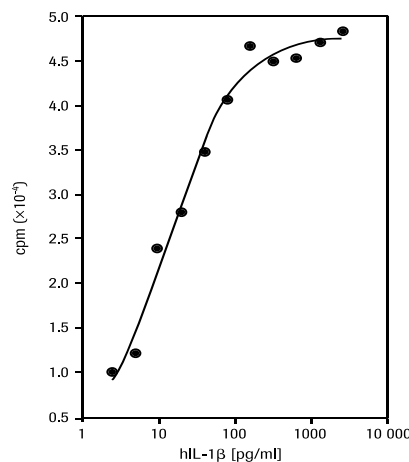


Fig. 1: [ $^3$ H]-thymidine incorporation into C3H/HeJ thymocytes in response to recombinant human interleukin-1 $\beta$  (hIL-1 $\beta$ ), using the method described.

## 4. Additional Information on this Product

### Background Information

Interleukin-1 (IL-1) is produced by a number of cell types, including activated macrophages, B-cells, fibroblasts, and keratinocytes (1,2). It mediates a wide range of biological activities, such as stimulation of thymocyte proliferation via induction of interleukin-2 (IL-2) release, stimulation of B-lymphocyte maturation and proliferation, fibroblast growth factor activity, and induction of acute-phase protein synthesis by hepatocytes. IL-1 has also been reported to stimulate prostaglandin and collagenase release from synovial cells, and to be identical to endogenous pyrogen and catabolin (1, 2-9). Two types of human interleukin-1 (hIL-1 $\alpha$  and hIL-1 $\beta$ ) have been described (1). Both types of hIL-1 stimulate proliferation and differentiation of T- and B-lymphocytes (7).

Recently, it has been shown that two different high-affinity IL-1 receptor molecules exist on different cell types. In both, human and mouse, T-cells and fibroblasts express an  $\approx$ 80 kDa receptor molecule, whereas an  $\approx$ 60 kDa receptor molecule was found on B-cells and neutrophils. The action of hIL-1 $\alpha$  and hIL-1 $\beta$  is mediated by both receptor molecules (8-11).

IL-1 appears to have a wide range of stimulatory effects on the maturation, differentiation, and growth of many cell types involved in inflammation and development. Cells whose growth is directly or indirectly stimulated by IL-1 are fibroblasts, synovial cells, endothelial cells, epithelial cells, bone marrow cells, T-lymphocytes, and B-lymphocytes (1, 2-6).

### Preparation

Recombinant Interleukin-1 $\beta$ , human (hIL-1 $\beta$ ) is produced in *E. coli* and purified by standard chromatographic techniques.

## Primary Structure

One polypeptide chain (154 amino acids), identical to natural human IL-1 $\beta$  (153 amino acids), but with an extra methionine at the amino-terminus (1,12,13).

## Purity

$\geq 95\%$  pure as determined by HPLC or SDS-PAGE [Endotoxin level:  $< 0.1$  EU/ $\mu$ g (LAL-test),  $< 10$  EU/ml (LAL-test)].

## Specific Activity

$> 5.0 \times 10^7$  U/mg; [ $^3$ H]-thymidine incorporation into mouse C3H/HeJ thymocytes in the presence of saturating amounts of hIL-2 (50 U/ml) (14-16) (hIL-1 $\beta$ , NIBSC 1<sup>st</sup> international standard, 86/680) (17) (see Figure 1).

## EC<sub>50</sub> Definition/Unit Definition

The amount of hIL-1b that is required to support half-maximal stimulation of DNA synthesis (BrdU incorporation) with mouse C3H/HeJ thymocytes in the presence of saturating amounts of hIL-2 (50 U/ml, 25 ng/ml) (1 unit equals  $< 0.02$  ng).

## Molecular Weight

17,000 Da

## Species Specificity

Recombinant IL-1 $\beta$ , human is effective on mouse and human cells.

## References

- 1 March, C. J. et al. (1985) *Nature* **315**, 641-647.
- 2 Dinarello, C.A. (1985) *J. Clin. Immunol.* **5**, 287-297.
- 3 Oppenheim, J.J. et al. (1986) *Immunol. Today* **7**, 45-56.
- 4 Durum, S.K., Schmidt, J.A. & Oppenheim, J.J. (1985) *Ann. Rev. Immunol.* **3**, 263-268.
- 5 Krakauer, T. (1986) *CRC Critical Rev. Immunol.* **6**, 213-244.
- 6 Nathan, C.F. (1987) *J. Clin. Invest.* **79**, 319-326.
- 7 Allison, A.C. (1986) *BioEssays* **3**, 260-263.
- 8 Matsushima, K. et al. (1986) *J. Immunol.* **136**, 4496-4501.
- 9 Chin, J. et al. (1987) *J. Exp. Med.* **165**, 70-86.
- 10 Chizzonite, R. (1989) *Proc. Natl. Acad. Sci. USA* **86**, 8029-8033.
- 11 Bomsztyk, K. et al. (1989) *Proc. Natl. Acad. Sci. USA* **86**, 8034-8038.
- 12 Wingfield, P. et al. (1986) *Eur. J. Biochem.* **160**, 491-497. Wingfield, P. et al. (1986) *Eur. J. Biochem.* **160**, 491-497.
- 13 Furutani, Y. et al. (1986) *Nucleic Acids Res.* **14**, 3167-3179.
- 14 Symons, J.A. et al. (1987) In: *Lymphokines and Interferons* (Clemens, M. J.; Morris, A.G. & Gearing, A.J.H., eds.) IRL Press, Oxford, Washington DC, pp. 269-289.
- 15 Remvig, L. et al. (1991) *Allergy* **46**, 59-67.
- 16 Falk, W.; Krammer, P.H. & Männel, D.N. (1987) *J. Immunol. Methods* **99**, 47-52.
- 17 Poole, S. & Gaines Das, R.E. (1991) *J. Immunol. Methods* **142**, 1-13.

\* available from Roche Applied Science

## 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 <b>1</b> , <b>2</b> , 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. |

### Ordering Information

Roche Applied Science offers a large selection of reagents and systems for life science research. For a complete overview of related products and manuals, please visit and bookmark our home page, [www.roche-applied-science.com](http://www.roche-applied-science.com).

### Disclaimer of License

Licensed from Cistron Biotechnology Inc. under U.S. Patents 4,766,069, 5,122,459 and 5,484,887.

### Regulatory Disclaimer

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

## Contact and Support

To ask questions, solve problems, suggest enhancements or report new applications, please visit our **Online Technical Support Site** at:

**[www.roche-applied-science.com/support](http://www.roche-applied-science.com/support)**

To call, write, fax, or email us, visit the Roche Applied Science home page, [www.roche-applied-science.com](http://www.roche-applied-science.com), and select your home country. Country-specific contact information will be displayed. Use the Product Search function to find Pack Inserts and Material Safety Data Sheets.



Roche Diagnostics GmbH  
Roche Applied Science  
68298 Mannheim  
Germany