

IL-1 beta

Interleukin-1 beta

human, recombinant, *E. coli*

Cat. No.	Amount
PR-401	10 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in PBS containing 50 μ g BSA/ μ g IL-1 β .

Purity

\geq 95% by SDS-PAGE

Description

Interleukins-1 (IL-1) α and β belong to the family of regulatory and inflammatory cytokines which play an important role in the regulation of acute inflammation. IL-1 β is involved inflammatory and immune responses and promotes wound healing. The closely related proteins IL-1 α and IL-1 β share approximately 62% homology in amino acid sequence but exhibit nearly identical biochemical action.

Selected References:

Bazan *et al.* (1996) A newly defined Interleukin-1. *Nature* **379**:591.
Arend *et al.* (1998) Interleukin-1 receptor antagonist - role in biology. *Annu. Rev. Immunol.* **16**:27.
Dunn (2002) Mechanisms by which cytokines signal the brain. *Neurobiology of the Immune System* **52**:43.