# IL-1β (soluble)(human),(recombinant)

Citations: 3

View Online »

# **Ordering Information**

Order Online »

ALX-522-056-C010	10µg
ALX-522-056-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

**View Online »** 

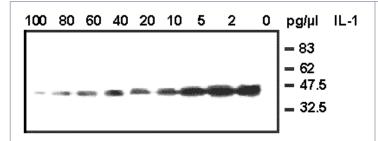


Figure: IL-1b activity was assessed by its ability to induce the activation of the NF-kB pathway. IkBa degradation was detected by Western blotting.

Method: Hela cells were grown to confluence in 6 well plates. IL-1b was diluted at the indicated concentrations in 500μl DMEM/Nut.Mix.F-12 10% fetal calf serum and added to the wells after being washed once with the same media. Cells were collected after 30 min and lysed. 100μg of cellular extract were separated on a SDS-PAGE and transferred on a nitrocellulose membrane. IkBa degradation was detected with a rabbit polyclonal antibody to IkBa and a goat anti-rabbit HRP-conjugated secondary antibody. Degradation of IkBa was detected at 10pg/μl IL-1b, under these conditions.

## **Handling & Storage**

**Use/Stability** Stable for at least 6 months after receipt when stored at -20°C.

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name IL-1F2, Interleukin-1β

**Concentration** 0.1mg/ml after reconstitution.

Endotoxin Content <0.1EU/µg purified protein (LAL test; Associates of CAPE

COD Inc.).

**Formulation** Lyophilized. Contains PBS.

Purity ≥90% (SDS-PAGE)

Reconstitution Reconstitute with 100µl sterile water.

**Source** Produced in *E. coli*. Human IL-1β (interleukin-1β) (aa 117-

270) is fused to a linker peptide (10 aa) and an N-terminal

FLAG<sup>®</sup>-tag.

**Specificity** Binds to human IL-1R (interleukin-1 receptor).

**Technical Info / Product Notes**Historical data has shown that IL-1β activates the NF-κB

pathway.

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID P01584



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com