

# IL-1 $\beta$ (soluble) (human), (recombinant)

Citations: 3

[View Online »](#)

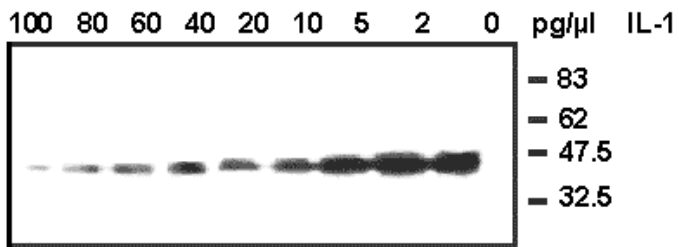
## Ordering Information

[Order Online »](#)

|                  |                          |
|------------------|--------------------------|
| ALX-522-056-C010 | 10 $\mu$ g               |
| ALX-522-056-3010 | SuperPack - 3x10 $\mu$ g |

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** IL-1b activity was assessed by its ability to induce the activation of the NF-κB pathway. IκBα 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. IκBα degradation was detected with a rabbit polyclonal antibody to IκBα and a goat anti-rabbit HRP-conjugated secondary antibody. Degradation of IκBα 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-1 $\beta$  |
| Concentration                  | 0.1mg/ml after reconstitution.   |
| Endotoxin Content              | <0.1EU/ $\mu$ g purified protein (LAL test; Associates of CAPE COD Inc.).  |
| Formulation                    | Lyophilized. Contains PBS.   |
| Purity                         | $\geq$ 90% (SDS-PAGE)  |
| Reconstitution                 | Reconstitute with 100 $\mu$ l sterile water.   |
| Source                         | Produced in <i>E. coli</i> . Human IL-1 $\beta$ (interleukin-1 $\beta$ ) (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-1 $\beta$ activates the NF- $\kappa$ B pathway.<br><br>FLAG is a registered trademark of Sigma-Aldrich Co.                         |
| UniProt ID                     | P01584   |



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)