

IL-1 receptor type I (human):Fc (human), (recombinant)

Ordering Information

[Order Online »](#)

ALX-522-057-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

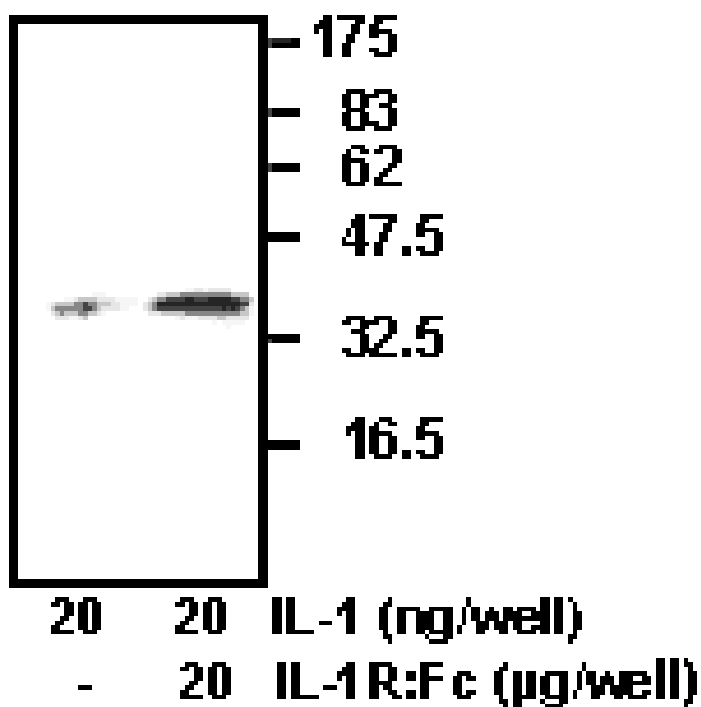


Figure: Inhibition of IL-1 (Prod. No. [ALX-522-056](#)) induced NF-κB activation by IL-1R:Fc. The degradation of IκBa upon IL-1 stimulation previously incubated with IL-1R:Fc was measured by Western blot. “”, IL-1 receptor type I (human):Fc (human)

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.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	IL-1RI, Interleukin-1 receptor type I, CD121a
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Lonza).
Formulation	Lyophilized. Contains PBS.
MW	~83kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 50μl sterile water. PBS containing at least 0.1% BSA should be used for further dilutions.
Source	Produced in HEK 293 cells. Human IL-1RI (interleukin-1 receptor type I) (aa 21-336) is fused to the Fc portion of human IgG1.
Specificity	Binds to human IL-1 (interleukin-1).
Technical Info / Product Notes	Historical data has shown that IL-1 receptor blocks IL-1 activity.
UniProt ID	P14778



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)