

IκBα (human), (recombinant) (untagged)

IκBα inhibits activation of the NF-κB pathway activation by associating with the major NF-κB heterodimeric complex p50/p65 and preventing its translocation from the cytoplasm into the nucleus, where it acts a transcription factor for genes involved in inflammation, autoimmune response, cell proliferation, and apoptosis. Inhibition of NF-κB pathway activation by IκBα is controlled by its sequential serine-phosphorylation, ubiquitination and proteasomal degradation. This process itself can be inhibited by SUMOylation of IκBα, providing an additional regulatory step for the activation of the NF-κB pathway.

Citations: 1

[View Online »](#)

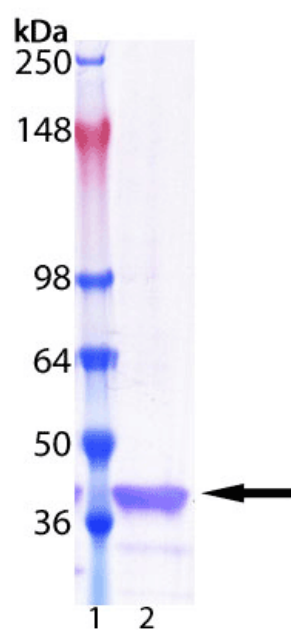
Ordering Information

[Order Online »](#)

BML-UW9975-0050	50μg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of Prod. No. BML-UW9975: Lane 1: MW marker, Lane 2: 1 μ g I κ B α (human), (recombinant) (untagged).

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

Uses:

1. NF-κB pathway studies, inhibition of NF-κB p50/p65, substrate for IκB kinase (IKK).
2. Investigation of IκBα phosphorylation/ubiquitinylation/SUMOylation.
3. Ubiquitin-proteasome mediated degradation studies.

Formulation

Liquid. In 50mM HEPES, pH 7.5, containing 200mM sodium chloride and 1mM DTT.

Source

Produced in *E. coli*.

UniProt ID

P25963



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)