

NF- κ B p65 subunit (human) (12-317), (recombinant) (untagged)

The eukaryotic nuclear factor κ B (NF- κ B) plays an important role in inflammation, autoimmune response, cell proliferation, and apoptosis by regulating the expression of genes involved in these processes.

Misregulation of the NF- κ B signal transduction pathway is observed in a variety of human cancers, especially those of lymphoid cell origin. The prototypical inducible NF- κ B is a heterodimeric complex consisting of p50 and p65, which normally exists in the cytoplasm of the cell as an inactive form due to association with its major inhibitory protein, I κ B α . I κ B α prevents the transport of NF- κ B into the nucleus by masking its nuclear localization signal. Activation of NF- κ B by extracellular inducers depends on the phosphorylation and subsequent degradation of I κ B α by the ubiquitin-proteasome pathway.

p65 subunit

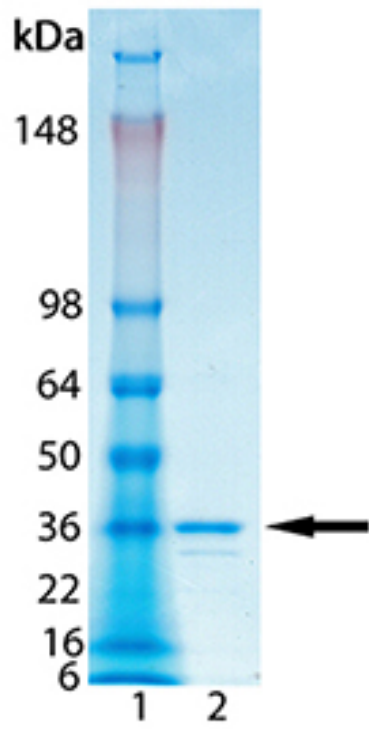
The NF- κ B p65 subunit is usually thought of as conferring the activation function of the NF- κ B complex, though its capacity to associate with κ B DNA in the absence of p50 and its potential utility as an independent transcription factor have also been demonstrated.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0 μg ; of purified NF- κB p65 Subunit (12-317) (human) (recombinant) (untagged) Protein Prod. No. BML-UW9995.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes	For NF-κB pathway studies.
Formulation	Liquid. In 1X Dulbecco's Phosphate Buffered Saline (DPBS), pH 7.4, containing 5% glycerol.
MW	~35kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> BL21 (λDE3). Recombinant human NF-κB p65 subunit (12-317).
UniProt ID	Q04206

Last modified: May 29, 2024



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)