

NF-κB p65 subunit (human) (12-317), (recombinant) (GST- tag)

The eukaryotic nuclear factor κB (NF-κB) 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-κB signal transduction pathway is observed in a variety of human cancers, especially those of lymphoid cell origin. The prototypical inducible NF-κB 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, IκBα. IκBα prevents the transport of NF-κB into the nucleus by masking its nuclear localization signal. Activation of NF-κB by extracellular inducers depends on the phosphorylation and subsequent degradation of IκBα by the ubiquitin-proteasome pathway.

p65 subunit

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

Citations: 2

[View Online »](#)

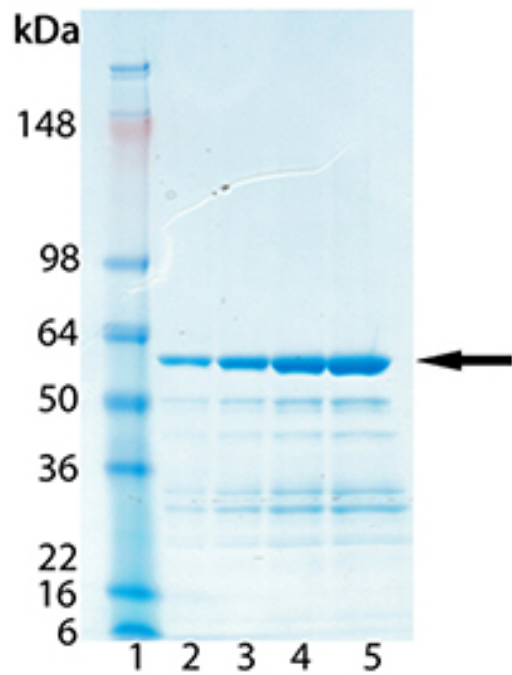
Ordering Information

[Order Online »](#)

BML-UW9990-0050	50μg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 0.5 μ g; Lane 3: 1.0 μ g; Lane 4: 2.0 μ g; Lane 5: 5.0 μ g of purified NF- κ B p65 Subunit (12-317) (human) (recombinant) (GST-tag) Protein Prod. No. BML-UW9990.

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 50mM HEPES, pH 7.5, containing 50mM sodium chloride and 1mM DTT.
MW	~60kDa
Purity	≥85% (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) is fused to a GST-tag.
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)