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

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, $I\kappa B\alpha$. $I\kappa B\alpha$ 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 $I\kappa B\alpha$ 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 »

Ordering Information

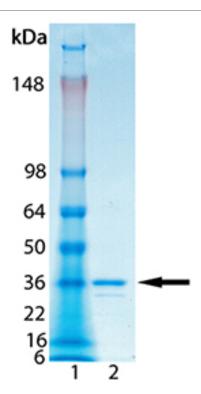
Order Online »

BML-UW9995-0050

50µg

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-kB 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 *E. coli* BL21 (λDE3). Recombinant human NF-

κB p65 subunit (12-317).

UniProt ID Q04206

Last modified: May 29, 2024

