

Recombinant NFκB1 p105 protein

Catalog No: 81143, 81843

Lot No: 10318001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.8 µg/µl

Source: Human

Buffer Contents: Recombinant NFκB1 / NFκB p105 protein is supplied in 25 mM HEPESNaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100 and 0.5 mM TCEP.

Background: NFκB1 (Nuclear Factor Kappa B Subunit 1), also known as NFKB1 or NF-kappa-B, is a pleiotropic transcription factor present in almost all cell types and is the endpoint of a series of signal transduction events that are initiated by a vast array of stimuli related to many biological processes such as inflammation, immunity, differentiation, cell growth, tumorigenesis and apoptosis. The 105 kD protein can undergo cotranslational processing by the 26S proteasome to produce a 50 kD protein. It is a Rel protein-specific transcription inhibitor and the 50 kD protein is a DNA binding subunit

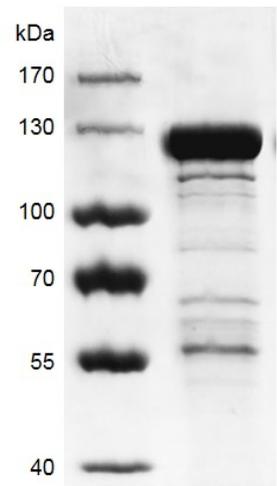
of the NF-kappa-B (NFKB) protein complex. NFKB1 is a homo- or heterodimeric complex formed by the Rel-like domaincontaining proteins RELA/p65, RELB, NFKB1/p105, NFKB1/p50, REL and NFKB2/p52 and the heterodimeric p65-p50 complex appears to be most abundant one. NFKB1 appears to have dual functions such as cytoplasmic retention of attached NF-kappa-B proteins by p105 and generation of p50 by a cotranslational processing. The proteasome-mediated process ensures the production of both p50 and p105 and preserves their independent function, although processing of NFκB1/p105 also appears to occur post-translationally.

Protein Details: Recombinant NF:kappa;B1 / NFκB p105 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_003989.2) with an N-terminal FLAG tag. The molecular weight of the protein is 107.1 kDa.

Application Notes: This protein is suitable for use in protein-protein interaction, in vitro transcription assay, binding assay.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

NFKB1 / NFκB p105



Recombinant NFKB1 / NFκB p105 protein gel

7.5% SDS-PAGE Coomassie staining
MW: 107.1 kDa
Purity: >85%