

NFKB2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # N13-31G

Lot # W308-2

Product Description

Recombinant human NFKB2 (1-454) was expressed in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_002502](#).

Gene Aliases

NFKB2-p52, NFKB2-p100, LYT10, LYT-10

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

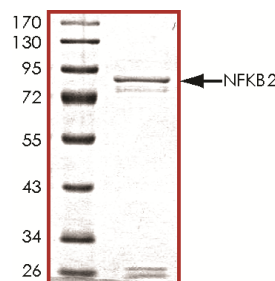
NFKB2 is a part of the NFKB complex that is present in various cell types that express cytokines, chemokines, growth factors, cell adhesion molecules, and some acute phase proteins in health and disease states. NFKB protein is activated by a wide variety of stimuli such as cytokines, oxidant-free radicals, inhaled particles, ultraviolet irradiation, and bacterial or viral products (1). Inappropriate activation of NFKB has been linked to inflammatory events associated with autoimmune arthritis, asthma, septic shock, lung fibrosis, glomerulonephritis, atherosclerosis, and AIDS (2). Overexpression of NFKB2 protects androgen sensitive LNCaP cells from apoptotic cell death and cell cycle arrest induced by androgen-deprivation.

References

1. Zhu, M. et al: NF-kappa-B2 is required for the establishment of central tolerance through an Aire-dependent pathway. *J. Clin. Invest.* 116: 2964-2971, 2006.
2. Tucker, E. et al: A novel mutation in the Nfkb2 gene generates an NF-kappa-B2 'super repressor' *J. Immun.* 179: 7514-7522, 2007.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity of NFKB2 protein was determined to be **>80%** by densitometry.
Approx. MW **84kDa**.

NFKB2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	N13-31G
Specific Lot Number	W308-2
Purity	>80%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.