

IκB-α (Ab-42) Antibody [#B7116]

Catalog Number: B7116-1, B7116-2, B7116

Amount: 50μg/50μl, 100μg/100μl, 200μg/200μl

Swiss-Prot No. : P25963

All Names: I-kappa-B-alpha, IKBA, IkappaBalpha, MAD3, Major histocompatibility complex enhancer-binding protein MAD3, NF-kappaB inhibitor alpha, NFKBI, NFKBIA, RL/IF-1

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Storage/Stability: Store at -20°C/1 year

Immunogen: The antiserum was produced against synthesized non-phosphopeptide derived from human IκB-α around the phosphorylation site of tyrosine 42 (E-E-Y^P-E-Q).

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity: IκB-α (Ab-42) antibody detects endogenous levels of total IκB-α protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000

References: Béraud C, et al. (1999) Proc Natl Acad Sci U S A 96(2): 429-434.

Sundström S, et al. (2005) J Virol 79(4): 2230-2239.

Liu L, et al. (1998) Mol Cell Biol 18(7): 4221-4234.

Shrivastava A, et al. (1998) J Virol 72(12): 9722-9728.

Franzoso G, et al. (1997) Genes Dev 11(24): 3482-3496.