

CNOT2 (Ab-101) Antibody [#B8387]

Catalog Number:	B8387
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZN8
Other Names:	CCR4-associated factor 2; CCR4-NOT transcription complex subunit 2; CCR4-NOT transcription complex, subunit 2; CDC36; CNOT2; HSPC131; MSTP046; NOT2 (negative regulator of transcription 2, yeast) homolog; NOT2H
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human CNOT2 around the phosphorylation site of serine 101 (S-L-S ^P -Q-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CNOT2 (Ab-101) antibody detects endogenous levels of total CNOT2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Albert T.K., Nucleic Acids Res. 28:809-817(2000). Zhang Q.-H., Genome Res. 10:1546-1560(2000). Ota T., Nat. Genet. 36:40-45(2004).