

NFAT3 (Ab-676) Antibody [#B0521]

Catalog Number:	B0521
Concentration:	1mg/ml
Swiss-Prot No.:	Q14934
Other Names:	NF-AT3, NF-ATc4, NFAC4, NFATC4, NFC4, Nuclear factor of activated T-cells, cytoplasmic 4, T cell transcription factor NFAT3, Transcription complex subunit NF-ATc4
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 NFAT3 around the phosphorylation site of serine 676 (K-R-S ^P -P-T).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NFAT3 (Ab-676) antibody detects endogenous levels of total NFAT3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Hoey T., Immunity 2:461-472(1995). Crabtree G.R., Cell 96:611-614(1999). Olsen J.V., Cell 127:635-648(2006).