

p47 phox Antibody [#B1161]

Catalog Number:	B1161
Concentration:	1mg/ml
Swiss-Prot No.:	P14598
Other Names:	47 kDa autosomal chronic granulomatous disease protein, 47 kDa neutrophil oxidase factor, NCF-1, NCF-47K, NCF1, Neutrophil NADPH oxidase factor 1, Neutrophil cytosol factor 1, Neutrophil cytosolic factor 1, P47 phox, p47-phox
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 peptide derived from Internal of human p47 phox.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p47 phox antibody detects endogenous levels of total p47 phox protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Volpp B.D., Proc. Natl. Acad. Sci. U.S.A. 86:7195-7199(1989). Volpp B.D., Proc. Natl. Acad. Sci. U.S.A. 86:9563-9563(1989). Lomax K.J., Science 245:409-412(1989).