

TTF2 Antibody [#C19640]

Catalog Number:	C19640
Concentration:	1mg/ml
Swiss-Prot No.:	O00358
Other Names:	Interferon-related developmental regulator 1, Nerve growth factor-inducible protein PC4
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 TTF2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TTF2 antibody detects endogenous levels of total TTF2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:10000
References:	Buanne P., Genomics 51:233-242(1998). Hillier L.W., Nature 424:157-164(2003). The MGC Project Team; Genome Res. 14:2121-2127(2004).