

TF2A1 Antibody [#C10702]

Catalog Number:	C10702
Concentration:	1mg/ml
Swiss-Prot No.:	P52655
Other Names:	Transcription initiation factor IIA subunit 1, General transcription factor IIA subunit 1, TFIIA-42, TFIIAL, Transcription initiation factor IIA alpha chain (TFIIA p35 subunit), Transcription initiation factor IIA beta chain (TFIIA p19 subunit), GTF2A1, TF2A1
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 C-terminal of human TF2A1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TF2A1 antibody detects endogenous levels of total TF2A1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ma D., Genes Dev. 7:2246-2257(1993). Dejong J., Genes Dev. 7:2220-2234(1993). Heilig R., Nature 421:601-607(2003).