

## TF2H2 Antibody [#C11010]

<b>Catalog Number:</b>	C11010
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q13888
<b>Other Names:</b>	TFIIH basal transcription factor complex p44 subunit, Basic transcription factor 2 44 kDa subunit, BTF2-p44, General transcription factor IIH polypeptide 2, GTF2H2, BTF2P44
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from N-terminal of human TF2H2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TF2H2 antibody detects endogenous levels of total TF2H2 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Humbert S., EMBO J. 13:2393-2398(1994). The MGC Project Team, Genome Res. 14:2121-2127(2004). van der Steege G., Eur. J. Hum. Genet. 3:87-95(1995).