

Tryptophan Hydroxylase (Ab-260) Antibody [#B1012]

Catalog Number:	B1012
Concentration:	1mg/ml
Swiss-Prot No.:	P17752
Other Names:	EC 1.14.16.4; TPH; TRPH; tryptophan 5-hydroxylase 1; tryptophan 5-monoxygenase 1
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 Tryptophan Hydroxylase around the phosphorylation site of serine 260 (H-S-S ^P -D-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Tryptophan Hydroxylase (Ab-260) antibody detects endogenous levels of total Tryptophan Hydroxylase protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Boulalard S., Nucleic Acids Res. 18:4257-4257(1990). Tipper J.P., Arch. Biochem. Biophys. 315:445-453(1994). The MGC Project Team; Genome Res. 14:2121-2127(2004).