

## Tryptophan Hydroxylase (Ab-58) Antibody [#B1011]

<b>Catalog Number:</b>	B1011
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P17752
<b>Other Names:</b>	EC 1.14.16.4, TPH, TRPH, tryptophan 5-hydroxylase 1, tryptophan 5-monoxygenase 1
<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 non-phosphopeptide derived from human Tryptophan Hydroxylase around the phosphorylation site of serine 58 (R-N-S <sup>P</sup> -E-F).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Tryptophan Hydroxylase (Ab-58) antibody detects endogenous levels of total Tryptophan Hydroxylase protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Ji Ying Sze, Development, Aug 2002; 129: 3901. Jeremy Veenstra-VanderWeele, Mol. Interv., Mar 2003; 3: 72. Nurgul Carkaci-Salli, J. Biol. Chem., Sep 2006; 281: 28105 - 28112.