

Tyrosine Hydroxylase (Ab-40) Antibody [#B0039]

- Catalog Number:** B0039-1, B0039-2, B0039
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q2M3B4
- All Names:** EC 1.14.16.2, TH isoform 3, TH isoform a, TH-4, TY3H, TYH, Tyrosine 3-hydroxylase, Tyrosine 5-monoxygenase, tyrosine hydroxylase
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human Tyrosine Hydroxylase around the phosphorylation site of serine 40 (R-Q-S^P-L-I).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Tyrosine Hydroxylase (Ab-40) antibody detects endogenous levels of total Tyrosine Hydroxylase protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1000 IHC: 1:50~100
IF: 1:500~1:1000 ELISA: 1:20000
- References:** Jahan Ara, PNAS, Jun 1998; 95: 7659.
E Carafoli, Crit. Rev. Biochem. Mol. Biol., Apr 2001; 36: 107.
P. William Conrad, J. Biol. Chem., Nov 1999; 274: 33709.
Y. Katayama, J Appl Physiol, Nov 1994; 77: 2086.