

PRODUCT INFORMATION



Tyrosine Hydroxylase (Phospho-Ser⁴⁰) Polyclonal Antibody

Item No. 10009414

Overview and Properties

Contents:	This vial contains 100 µl of affinity-purified polyclonal antibody.
Synonym:	TH
Immunogen:	Synthetic peptide from amino acid residues surrounding phospho-Ser ⁴⁰ of rat TH
Molecular Weight:	60 kDa
Cross Reactivity:	(+) Mammalian TH, Non-mammalian TH
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
Host:	Rabbit
Isotype:	IgG
Applications:	Immunofluorescence (IF), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:1,000 for IF, IHC, and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

Tyrosine hydroxylase (TH) is the rate-limiting enzyme in the synthesis of the catecholamines dopamine and norepinephrine. TH antibodies can therefore be used as markers for dopaminergic and noradrenergic neurons in a variety of applications including depression, schizophrenia, Parkinson's disease, and drug abuse.¹⁻³ TH antibodies can also be used to explore basic mechanisms of dopamine and norepinephrine signaling.⁴⁻⁶ The activity of TH is also regulated by phosphorylation.⁷⁻⁹ Phospho-specific antibodies for the phosphorylation sites on TH can be used to great effect in studying this regulation and in identifying the cells in which TH phosphorylation occurs.

References

1. Kish, S.J., Kalasinsky, K.S., Derkach, P., *et al.* *Neuropsychopharmacology* **24(5)**, 561-567 (2001).
2. Zhu, M.-Y., Klimek, V., Haycock, J.W., *et al.* *J. Neurosci. Methods* **99(1-2)**, 37-44 (2000).
3. Zhu, M.-Y., Klimek, V., Dilley, G.E., *et al.* *Biol. Psychiatry* **46(9)**, 1275-1286 (1999).
4. Witkovsky, P., Gabriel, R., Haycock, J.W., *et al.* *J. Chem. Neuroanat.* **19(2)**, 105-116 (2000).
5. Salvatore, M.F., Waymire, J.C., and Haycock, J.W. *J. Neurochem.* **79(2)**, 349-360 (2001).
6. Dunkley, P.R., Bobrovskaya, L., Graham, M.E., *et al.* *J. Neurochem.* **91(5)**, 1025-1043 (2004).
7. Haycock, J.W., Ahn, N.G., Cobb, M.H., *et al.* *Proc. Natl. Acad. Sci. USA* **89(6)**, 2365-2369 (1992).
8. Haycock, J.W., Bennett, W.F., George, R.J., *et al.* *J. Biol. Chem.* **257(22)**, 13699-13703 (1982).
9. Jedynak, J.P., Ali, S.F., Haycock, J.W., *et al.* *J. Neurochem.* **82(2)**, 382-388 (2002).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/29/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM