

# Product Information



## Dopamine $\beta$ -hydroxylase (C-Term; human) Polyclonal Antibody

Item No. 10009370

<b>Synonym:</b>	DBH
<b>Supplied as:</b>	100 $\mu$ l of affinity-purified antibody in 10 mM (pH 7.5), 150mM NaCl, 100 $\mu$ g per ml BSA and 50% glycerol.
<b>Host:</b>	Sheep
<b>Antigen:</b>	Peptide from the C-terminal region of human DBH
<b>Cross Reactivity:</b>	(+) Human, mouse, and non-human primate DBH
<b>Stability:</b>	$\geq$ 1 year at $-20^{\circ}\text{C}$
<b>Application:</b>	The recommended starting dilution for western blot is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Dopamine  $\beta$ -hydroxylase (DBH) catalyzes the conversion of dopamine to norepinephrine and serves as a marker of noradrenergic cells. DBH antibodies and antibodies for other markers of catecholamine biosynthesis are widely used as markers for dopaminergic and noradrenergic neurons in a variety of applications including depression, schizophrenia, Parkinson's disease, and drug abuse.<sup>1-3</sup> The expression of DBH is also elevated during stress.<sup>4</sup>

### References

1. Kish, S.J., Kalasinsky, K.S., Derkach, P., *et al.* Striatal dopaminergic and serotonergic markers in human heroin users. *Neuropsychopharmacology* **24**, 561-567 (2001).
2. Zhu, M.Y., Klimek, V., Haycock, J.W., *et al.* Quantitation of tyrosine hydroxylase protein in the locus coeruleus from postmortem human brain. *J. Neurosci. Meth.* **99**, 37-44 (2000).
3. Zhu, M.Y., Klimek, V., Diley, G.E., *et al.* Elevated levels of tyrosine hydroxylase in the locus coeruleus in major depression. *Biol. Psychiatry* **46**, 1275-1286 (1999).
4. Sabban, E.L. and Kvetnansky, R. Stress-triggered activation of gene expression in catecholaminergic systems: Dynamics of transcriptional events. *Trends Neurosci.* **24**, 91-98 (2001).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10009370](http://www.caymanchem.com/catalog/10009370)

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 12/16/2011