

Product Information



Dopamine Transporter (C-Term) Polyclonal Antibody

Item No. 10009372

Synonym:	DAT (C-Term)
Supplied as:	100 µl of peptide affinity-purified antibody in 10 mM HEPES, pH 7.5, 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.
Host:	Rabbit
Antigen:	Peptide from the C-terminal region of human DAT
Cross Reactivity:	Western blot (WB): (+) human and mouse striatal samples (SDS-solubilized) DAT; Immunohistochemistry (IHC): (+) Macaque monkey brain DAT
Stability:	≥1 year at -20°C
Applications:	The recommended starting dilution for WB and IHC is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

The dopamine transporter (DAT) is responsible for the reaccumulation of dopamine after it has been released. DAT 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.¹⁻³ Levels of DAT protein expression are altered by chronic drug administration.⁴

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. Wilson, J.M., Kalasinsky, K.S., Levey, A.I., *et al.* Striatal dopamine nerve terminal markers in human, chronic methamphetamine users. *Nat. Med.* **2**, 699-703 (1996).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009372

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

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

Web

www.caymanchem.com