

# Product Information



## HDAC11 Polyclonal Antibody

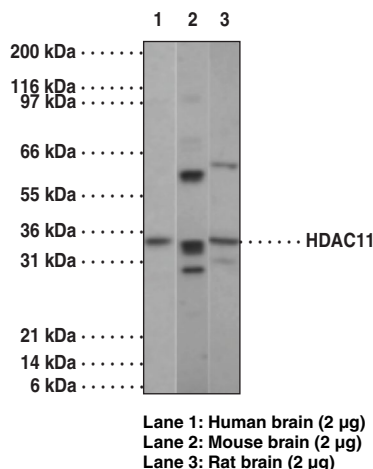
Item No. 13504

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.02% gelatin and 0.05% sodium azide.
- Synonym:** Histone Deacetylase 11
- Antigen:** Synthetic peptide from human HDAC11 amino acids 182-199
- Host:** Rabbit
- Cross Reactivity:** (+) Human, mouse, and rat HDAC11
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** Western Blot: 1-3 µg/ml

Histone deacetylase (HDAC) and histone acetyltransferase (HAT) are enzymes that regulate transcription by selectively deacetylating or acetylating the ε-amino groups of lysines located near the amino termini of core histone proteins.<sup>1</sup> Eleven members of HDAC family have been identified in the past several years.<sup>2,3</sup> These HDAC family members are divided into two classes, I and II. The newest member of this family, HDAC11, has been cloned by Gao *et al.*<sup>4</sup> HDAC11 transcripts were limited to kidney, heart, brain, skeletal muscle, and testis. These results suggested that each member of the HDAC family exhibits a different, individual substrate specificity and function *in vivo*.

### References

1. Cress, W.D. and Seto, E.J. Histone deacetylases, transcriptional control, and cancer. *J. Cell. Physiol.* **184**, 1-16 (2000).
2. Meinke, P.T., Liberator, P. Histone deacetylase: A target for antiproliferative and antiprotozoal agents. *Curr. Med. Chem.* **8(2)**, 211-235 (2001).
3. Nakayama, T., Takami, Y. Participation of histones and histone-modifying enzymes in cell functions through alterations in chromatin structure. *J. Biochem.* **129**, 491-499 (2001).
4. Gao, L., Cueto, M.A., Asselbers, F., *et al.* Cloning and functional characterization of HDAC11, a novel member of the human histone deacetylase family. *J. Biol. Chem.* **277(28)**, 25748-25755 (2002).



### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13504](http://www.caymanchem.com/catalog/13504)

### 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](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, 01/21/2013