

Product Information

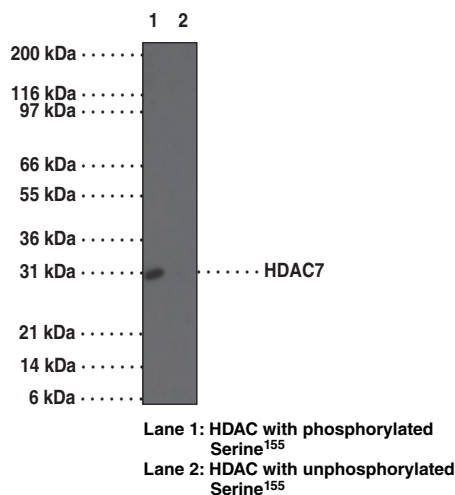


HDAC7 (Phospho-Ser¹⁵⁵) Polyclonal Antibody

Item No. 13500

- Contents:** This vial contains 50 µg of peptide-purified IgG in 100 µl of PBS containing 0.05% BSA and 0.05% sodium azide.
- Synonym:** Histone Deacetylase 7
- Antigen:** Synthetic peptide from human HDAC7 containing phospho-Ser¹⁵⁵
- Host:** Rabbit
- Cross Reactivity:** (+) Chimpanzee, bovine, canine, human, monkey, mouse, and rat HDAC7
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 0.1-1 µg/ml.

Histone deacetylase 7 (HDAC7) is a thymus-specific, intracellular transcription factor belonging to the class IIa HDAC family. Reports suggest that dephosphorylation and nuclear localization of HDAC7 is promoted by myosin phosphatase, thus regulating the repression of Nur77 and inhibition of apoptosis in CD4⁺/CD8⁺ double-positive thymocytes. HDAC7 plays a specific role in maintaining vascular integrity by repressing the expression of matrix metalloproteinase (MMP) 10. It promotes repression mediated by transcriptional corepressor NCOR2 and is an efficient corepressor of the androgen receptor (AR). It is also responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3, and H4).



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13500

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, 01/29/2013

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