

Product Information

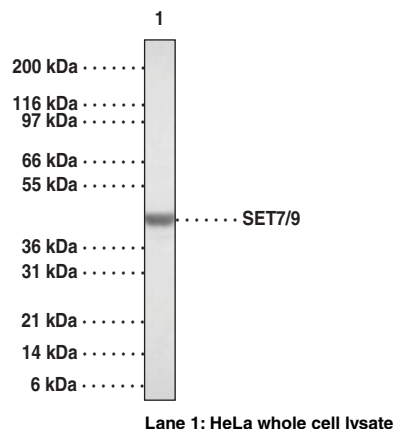


SET7/9 Polyclonal Antibody

Item No. 13731

Contents:	This vial contains 200 µl of protein G-purified IgG in 200 µl PBS containing 0.2% gelatin and 0.05% sodium azide.
Antigen:	Human SET7/9 amino acids 131-145 (GEVNEDGEMTGEKIA) and 336-352 (GYDHSPPGKSGPEAPEW)
Host:	Rabbit
Cross Reactivity:	(+) Human and mouse SET7/9
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Application:	Western blot (WB). The recommended starting dilution for WB is 1:500 - 1:1,000.

Diverse signal transduction pathways impinging on the N-terminal tails of histones lead to a number of post-translational modifications including acetylation, phosphorylation, poly(ADP-ribosylation), ubiquitination, and methylation. These modifications play critical roles in regulating chromatin structure and gene expression. Histone methyltransferases (HMTases) selectively methylate evolutionarily conserved arginine or lysine residues, primarily in the N-terminal tails of histones H3 and H4. SET7/9 is a histone specific HMTase that methylates histone H3 lysine 4. SET7/9 transfers methyl groups to lysine 4 of histone H3 in complex with S-adenosyl-L-methionine. Human SET7/9 is a 366 amino acid protein.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13731

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

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, 02/06/2013