

Product Information

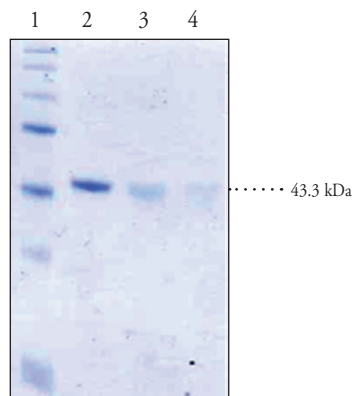
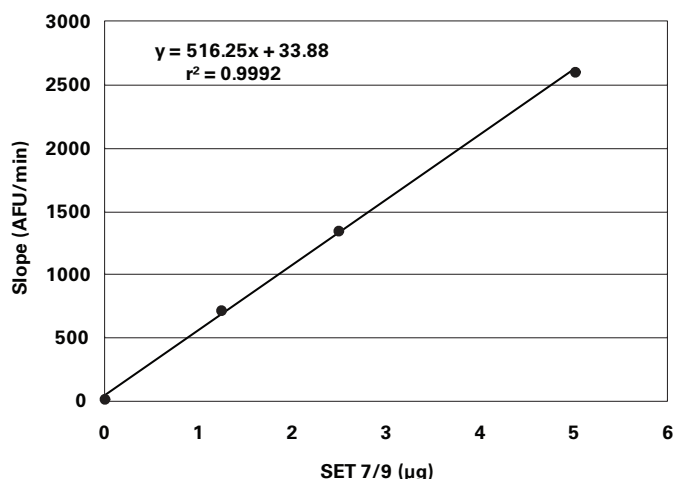


SET7/9 (human recombinant)

Item No. 10320 • Lot No. XXXX

Synonyms:	KMT7, SETD7/9, SET Domain-Containing Protein 7/9
Source:	Recombinant N-terminal His-tagged SET7/9 purified from <i>E. coli</i>
Amino Acids:	1-366 (Full-length)
M_r:	43.3 kDa
Purity:	≥95%
Stability:	≥6 months at -80°C
Supplied in:	50 mM sodium phosphate, pH 7.2, containing 100 mM sodium chloride and 20% glycerol
Activity:	XX U/ml. Determined using 50 μM human TAF10 peptide (amino acids 186-195) (Item No. 10228) (Ac)SKSKDRKYTL at 37°C using Cayman's Methyltransferase Fluorometric Assay Kit (Item No. 700150).
Specific Activity:	XX μmoles/min/mg
Protein Concentration:	XX mg/ml

Methylation of lysine can promote transcriptional activation or repression and is critical in regulating histone function.¹ Lysine residues can be mono-, di-, or tri-methylated, and unlike most SET proteins, SET7/9 is exclusively a mono-methylase.² SET7/9 methylates histone H3, tumor suppressor p53, and transcription factor TAF10.³ Recently, SET7/9 was shown to catalyze methylation of p53 in response to DNA damage thereby activating p53 for subsequent acetylation.¹ SET7/9 is able to modulate p53 activity in a human cancer cell line, implying that it may play a significant role in human tumorigenesis.



1. MW Markers
2. SET7/9 (4 μg)
3. SET7/9 (1 μg)
4. SET7/9 (0.5 μg)
Coomassie stained SDS-PAGE of Purified SET7/9.

References

1. Kurash, J.K., Lei, H., Shen, Q., *et al. Molecular Cell* **29**, 392-400 (2008).
2. Xiao, B., Jing, C., Wilson, J.R., *et al. Nature* **421**, 652-656 (2003).
3. Couture, J.-F., Collazo, E., Hauk, G., *et al. Nature Structural and Molecular Biology* **13(2)**, 140-146 (2006).

Related Products

TAF 10 Peptide - Item No. 10228 • JMJD2D (human recombinant) - Item No. 10335 • PRMT1 (human recombinant) - Item No. 10350 • G9a (human recombinant, partial) - Item No. 10353 • Methyltransferase Colorimetric Assay Kit - Item No. 700140 • Methyltransferase Fluorometric Assay Kit - Item No. 700150

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, 04/05/2011

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