

PRODUCT INFORMATION



Methylated Lysine Polyclonal Antibody HRP Conjugate

Item No. 13729

Overview and Properties

Contents:	This vial contains 400 µl of protein A-purified polyclonal antibody.
Immunogen:	Methylated KLH
Species Reactivity:	(+) Species independent
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol
Concentration:	0.25 mg/ml
Host:	Rabbit
Applications:	ELISA, Immunocytochemistry (ICC), Immunofluorescence (IF), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution for is 1:50 for ICC and IF and 1:5,000 for WB; other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

Post-translational modifications of proteins play critical roles in the regulation and function of many known biological processes. Proteins can be post-translationally modified in many different ways, and a common post-transcriptional modification of lysine involves methylation.¹ Lysine can be methylated once, twice, or three times by lysine methyltransferases. The transfer of methyl groups from S-adenosyl methionine to histones is catalyzed by enzymes known as histone methyltransferases. Histones which are methylated on certain residues can act epigenetically to repress or activate gene expression.^{1,2}

References

1. Yang, X.J. Multisite protein modification and intramolecular signaling. *Oncogene* **24(10)**, 1653-1662 (2005).
2. Melcher, M., Schmid, M., Aagaard, L., *et al.* Structure-function analysis of SUV39H1 reveals a dominant role in heterochromatin organization, chromosome segregation, and mitotic progression. *Mol. Cell Biol.* **20(10)**, 3728-3841 (2000).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/15/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM