

PRODUCT INFORMATION



Methylated Lysine Polyclonal Antibody

Item No. 31686

Overview and Properties

| | |
|--------------------------|--|
| Contents: | This vial contains 100 µg of protein A-affinity purified polyclonal antibody. |
| Immunogen: | Methylated KLH |
| Cross Reactivity: | (+) Mono- and di-methylated lysine residues in multiple species (cross-reactivity with tri-methyl lysine has not been tested); (-) Acetylated lysine |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | PBS with 50% glycerol |
| Concentration: | 1 mg/ml |
| Host: | Rabbit |
| Isotype: | IgG |
| Applications: | ELISA, immunohistochemistry (IHC), immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution for WB is (1:2,000-1:5,000). ELISA, IHC, IP and 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, 03/09/2023

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