

PRODUCT INFORMATION



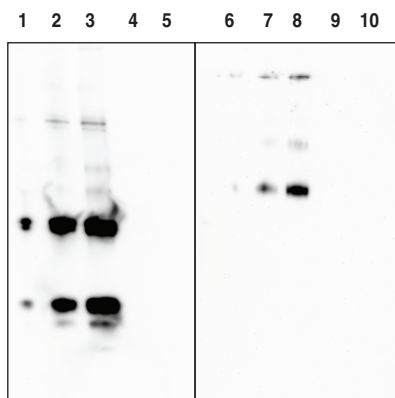
Diglycyl-Lysine Monoclonal Antibody (Clone GX41)

Item No. 21096

Overview and Properties

Contents:	This vial contains 100 µg of protein G-purified monoclonal antibody.
Immunogen:	Diglycine-modified histone peptide
Cross Reactivity:	(+) Diglycyl-modified lysine; (-) Unmodified lysine
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Clone:	GX41
Host:	Mouse
Isotype:	IgG1
Applications:	Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution for WB is 1:1,000. The recommended conditions for IP are 2-5 µg of antibody per test. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Calf Thymus Histone H2A-Gly-Gly (10 ng)
Lane 2: Calf Thymus Histone H2A-Gly-Gly (50 ng)
Lane 3: Calf Thymus Histone H2A-Gly-Gly (100 ng)
Lane 4: Calf Thymus Histone H2A (0.5 µg)
Lane 5: Calf Thymus Histone H2A (1 µg)
Lane 6: BSA-Gly-Gly (10 ng)
Lane 7: BSA-Gly-Gly (50 ng)
Lane 8: BSA-Gly-Gly (100 ng)
Lane 9: BSA (0.5 µg)
Lane 10: BSA (1 µg)

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, 08/16/2024

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

PRODUCT INFORMATION



Description

Diglycyl-lysine is a ubiquitin remnant-containing peptide.^{1,2} Diglycyl-lysine is formed during protein digestion with trypsin, which cleaves the arginine residue of ubiquitin, leaving diglycyl conjugated to lysine in protein digests.² It has been used as a marker of prior ubiquitin conjugation in trypsin protein digests.¹ Cayman's Diglycyl-Lysine Monoclonal Antibody (Clone GX41) can be used for Western blot, and immunoprecipitation applications.

References

1. Xu, G. and Jaffrey, S.R. Proteomic identification of protein ubiquitination events. *Biotechnol. Genet. Eng. Rev.* **29(1)**, 73-109 (2013).
2. Xu, G., Paige, J.S., and Jaffrey, S.R. Global analysis of lysine ubiquitination by ubiquitin remnant immunoaffinity profiling. *Nat. Biotechnol.* **28(8)**, 868-873 (2010).

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