

Diglycyl Lysine monoclonal antibody (GX41)

Highly specific monoclonal for ubiquitin remnant profiling via immunoprecipitation.

Protein ubiquitination is a fundamentally important post-translational modification that regulates myriad cell signaling pathways in eukaryotes. Alterations in protein ubiquitination have been linked to cancers and neurodegenerative, cardiovascular, immunological, and inflammatory diseases.

Identification of exact ubiquitylation sites is challenging under physiological conditions, but the recent development of an immunoprecipitation approach allows for enrichment and identification of peptides derived from the ubiquitinated portion of proteins.

When ubiquitinated protein is digested with trypsin a diglycine residue remains at any lysine site of ubiquitin attachment. Diglycyl Lysine, mAb (GX41) facilitates easy capture and enrichment of peptides derived from ubiquitinated proteins by recognizing the diglycyl-lysine residue with no detectable reactivity to non modified lysines.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-908-310-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store unopened vial at -20°C until required for use. Store diluted antibody at 2-4°C and use within 1 month. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IP, WB
Clone	GX41
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Mouse
Immunogen	Diglycine-modified histone from calf thymus, Type III-S, lysine rich fraction.
Isotype	IgG1κ
Purity Detail	Protein G purified.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Species independent
Specificity	Recognizes diglycyl modified lysine, a ubiquitin remnant detectable after trypsin digestion.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)