

V5-tag polyclonal antibody

Epitope tagging is a powerful and versatile strategy for detecting and purifying proteins expressed by cloned genes. To utilize this feature, protein expression vectors are typically engineered with a nucleotide sequence that encodes the epitope tag. The gene of interest is cloned in-frame relative to the tag and, upon expression, the protein of interest is synthesized as a fusion protein with the epitope tag. Fusion protein detection and/or purification is mediated by highly specific antibodies to the engineered protein, thus eliminating the need for antibodies to proteins from each newly cloned gene. Commonly used epitope tags include glutathione-S-transferase (GST), c-myc, 6-histidine (6X-His), FLAG, green fluorescent protein (GFP), maltose binding protein (MBP), influenza A virus haemagglutinin (HA), b-galactosidase, and GAL4.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to GKPIP NPLLGLDST.

Isotype IgG

Purity Detail Antigen affinity chromatography

Species Reactivity Species independent

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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