

Vesicular stomatitis virus (VSV-G) tag polyclonal antibody

Epitope tagging is a widely accepted technique that fuses an epitope peptide to a certain protein as a marker for gene expression. With this technique, gene expression can be easily monitored on Western blotting, immunoprecipitation and immunofluorescence utilizing an antibody that recognizes such an epitope. The VSV-G Tag (YTDIEMNRLGK) corresponds to the partial peptide sequence of the vesicular stomatitis virus glycoprotein.

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-MSA-115-E	100µl
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

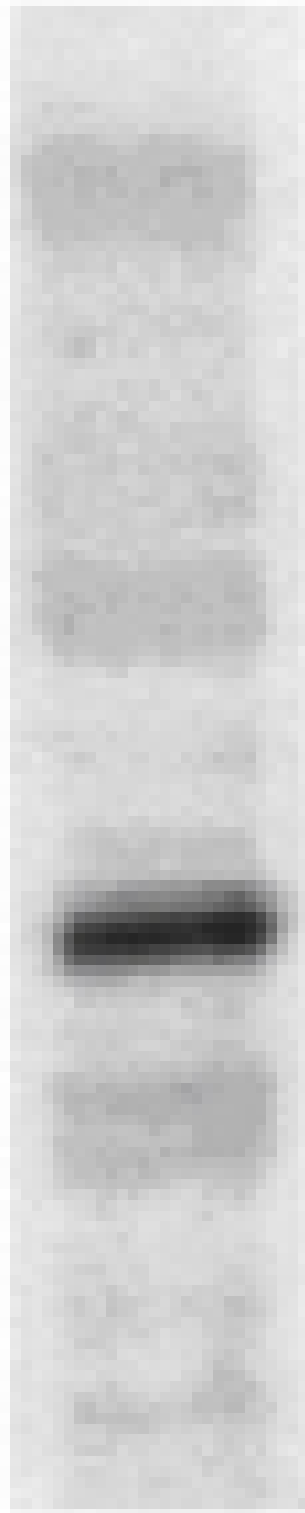
kDa

97 -

66 -

45 -

30 -



Western blot analysis of VSV-G-tag pAb expression in transfectant using 563

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, WB

Formulation Liquid. In PBS containing 50% glycerol, pH 7.2.

Host Rabbit

Immunogen Vesicular stomatitis virus (VSV-G) tag (YTDIEMNRLGK) coupled to GST.

Purity Detail Ammonium sulfate precipitation purified.

Recommendation Immunocytochemistry (10µg/ml)Western Blot (1:100, colorimetric)Suggested
Dilutions/Conditions dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

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



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)