

Rabies virus monoclonal antibody (Rab-50)

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

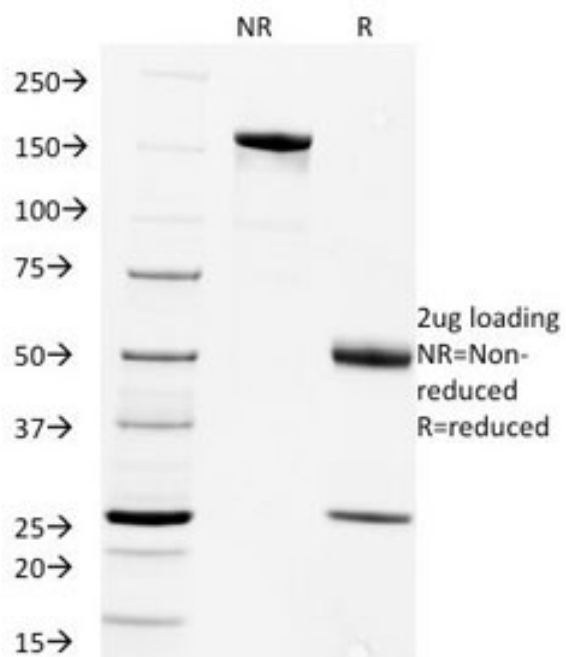
Ordering Information

[Order Online »](#)

ENZ-ABS914-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis of Purified, BSA-Free Rabies Virus Antibody (clone Rab-50). Confirmation of Integrity and Purity of the Antibody.

Handling & Storage

Handling	Store the Rabies Virus antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide).
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, ICC, IF
Clone	Rab-50
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Mouse
Immunogen	Purified virions of SAD-Vnukovo strain followed by Pitman-Moore strain.
Isotype	IgG2bk
Purity Detail	Protein G affinity chromatography purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (Acetone-fixed frozen) (0.5-1µg/ml for 30 min)Immunofluorescence (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Rabies Virus of strains SAD-Vnukovo and Pitman Moore

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Last modified: July 12, 2024
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-
be@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