

Rabies virus monoclonal antibody (Rab-50)

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

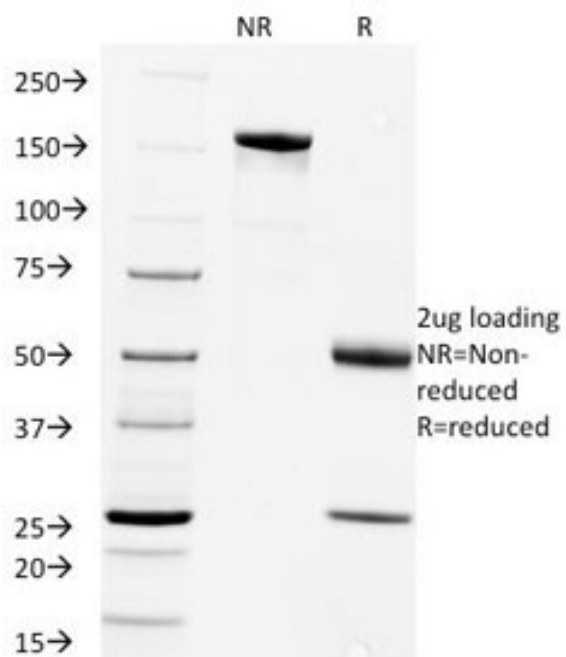
Ordering Information

[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.
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-be@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)