

PMEL17 recombinant monoclonal antibody (PMEL/1825R)

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

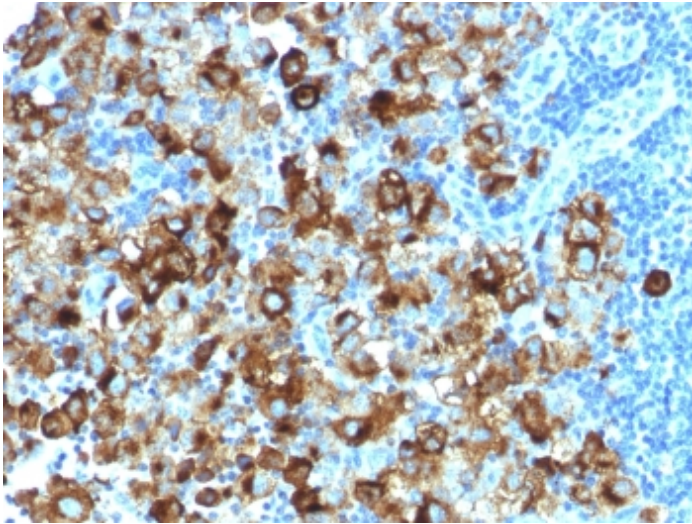
Ordering Information

[Order Online »](#)

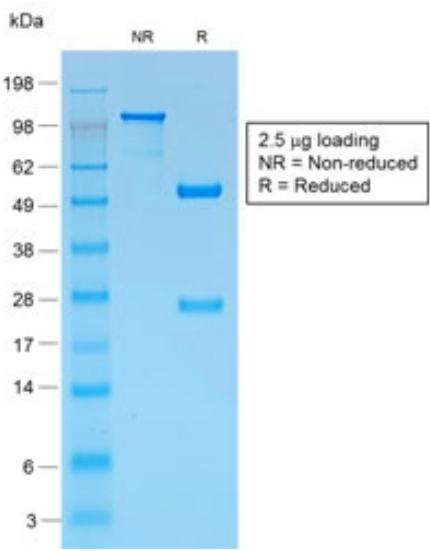
ENZ-ABS911-0100	100µg
-----------------	-------

Manuals, SDS & CofA

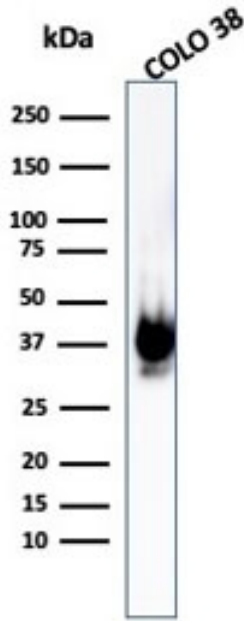
[View Online »](#)



IHC testing of FFPE human melanoma with recombinant gp100 antibody (clone PMEL/1825R). Required HIER: boil tissue sections in 10mM citrate buffer, pH 6, for 10-20 min followed by cooling at RT for 20 min.



SDS-PAGE analysis of purified, BSA-free recombinant gp100 antibody (clone PMEL/1825R) as confirmation of integrity and purity.



Western blot testing of human COLO-38 cell lysate with recombinant gp100 antibody (clone PMEL/1825R).

Handling & Storage

Handling	Store the recombinant gp100 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

Alternative Name	Melanoma lineage-specific antigen gp100, Melanoma-associated ME20 antigen
Application	IHC (PS), WB
Clone	PMEL/1825R
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Recombinant human gp100.
Isotype	IgGk
Purity Detail	Protein A affinity chromatography purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (FFPE) (0.5-1µg/ml for 30 min at RT)Western blot (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P40967

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)