

Melanoma marker (human) monoclonal antibody (HMB45)

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

Citations: 22

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-C34930	~10 µg - 0.5ml
------------	----------------

Manuals, SDS & CofA

[View Online »](#)

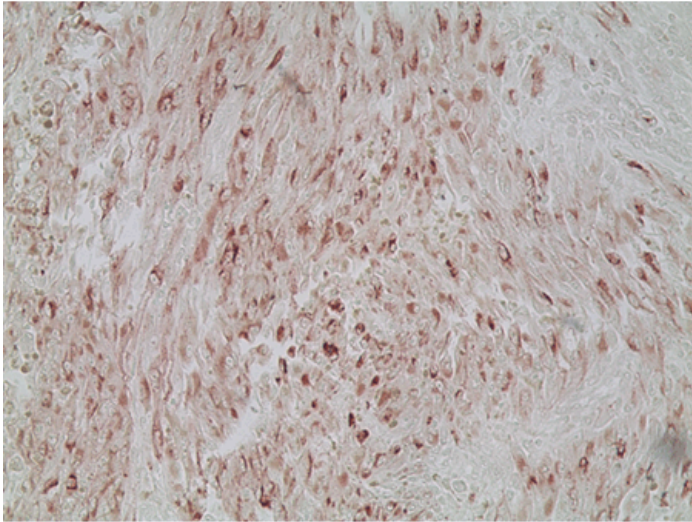


Figure 2: Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45), (Prod. No. ENZ-C34930) at a dilution of 1:20.

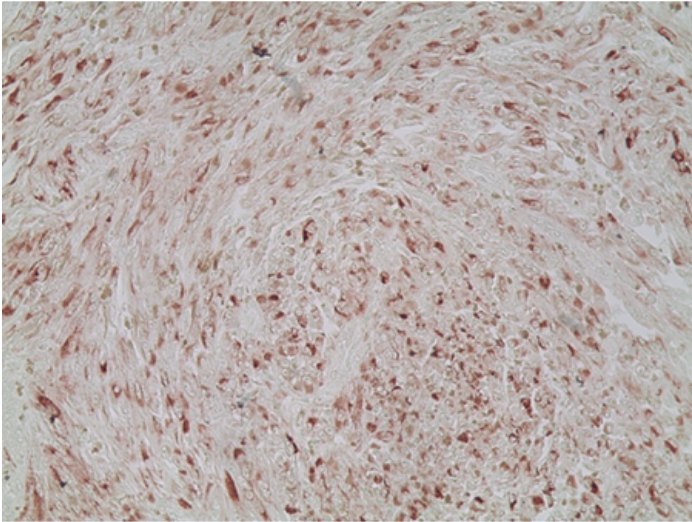


Figure 1: Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45), (Prod. No. ENZ-C34930) at a dilution of 1:20.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IHC (PS)
Clone	HMB45
Crossreactivity	Does not cross-react with normal melanocytes, intradermal nevus cells or other neoplasms including sarcomas, carcinomas and glial tumors.
Formulation	Liquid. In PBS containing 0.1% sodium azide and 1% BSA.
Host	Mouse
Immunogen	Human axillary lymph node tissue containing pigmented melanoma metastases.
Isotype	IgG1
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections; this antibody can be used for staining of paraffin-embedded tissue sections fixed in neutral buffered formalin and is compatible with most commonly used histological fixatives. Choose a good quality enzyme-linked anti-mouse IgG1 antibody for detection). A 20- to 50-fold dilution of the antibody is suggested. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually.
Source	Purified from ascites.
Species Reactivity	Human

Specificity

Recognizes a human melanoma-specific antigen present in melanocytic cells and junctional nevi. Does not cross-react with normal melanocytes, intradermal nevus cells or other neoplasms including sarcomas, carcinomas and glial tumors.

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)