

Melanoma marker (human) monoclonal antibody (HMB45)

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

Citations: 23

[View Online »](#)

Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)

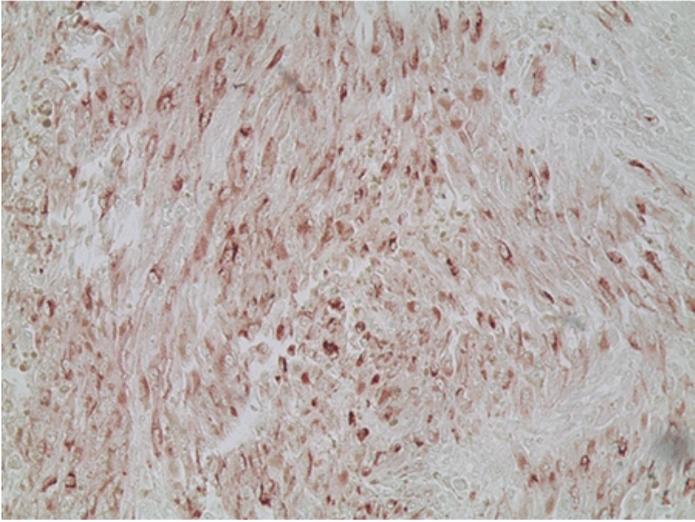


Figure 2: Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45), (Prod. No. ENZ-C34930) at a dilution of 1:20.

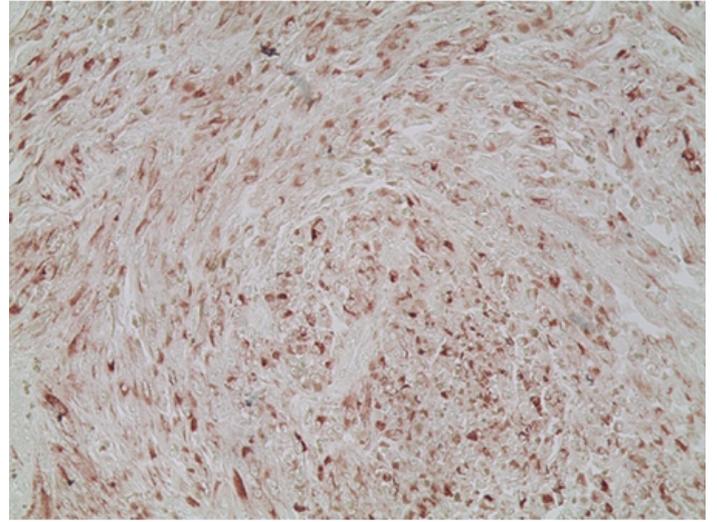


Figure 1: Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45), (Prod. No. ENZ-C34930) at a dilution of 1:20.

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](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
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