

Melanoma marker (human) monoclonal antibody (HMB45) (Ready-to-Use)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|-----------|-------------|
| ENZ-30930 | ~6 µg - 6ml |
|-----------|-------------|

Manuals, SDS & CofA

[View Online »](#)

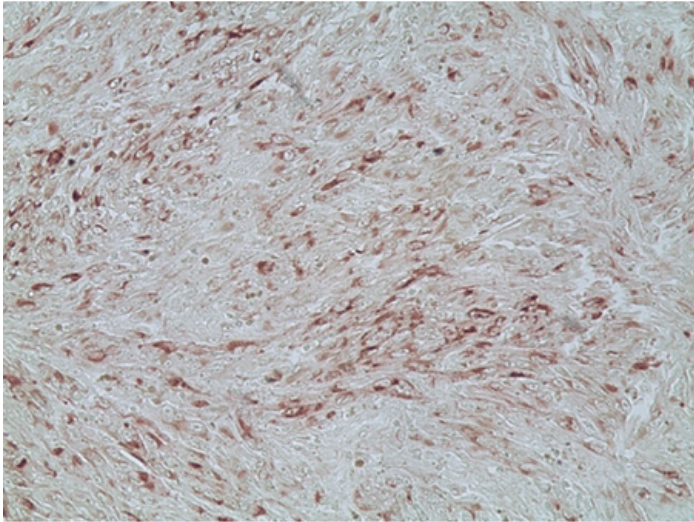


Figure 1. Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45) (Ready-To-Use), (Prod. No. ENZ-30930). Antibody is applied as is.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°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. *Ready-to-use* solution 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). 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



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-fr@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

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee**.