

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 »](#)

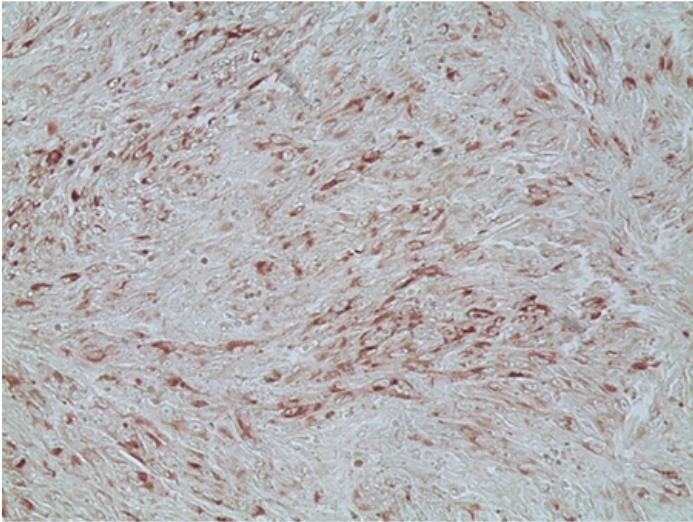


Figure 1. Immunohistochemistry analysis of human melanoma skin tissue using Melanoma marker (human) mAb (HMB45) (Ready-To-Use), (Prod. No. ENZ-30930). Antibody is applied as is.

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. <i>Ready-to-use</i> 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.
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
---	--	--	--	---	---