

Nucleolin monoclonal antibody (4E2)

Nucleolin (identical to human DNA helicase IV) is an abundant, non-ribosomal protein of the nucleolus involved in the organization of nuclear chromatin, rDNA transcription, packaging of pre-RNA, ribosome assembly, nucleocytoplasmic transport, cytokinesis, nucleogenesis, and apoptosis. A predominantly nuclear protein, the presence of nucleolin at the surface of cancer cells suggests that it could be valuable as a marker for the diagnosis of cancer.

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

Citations: 5

[View Online »](#)

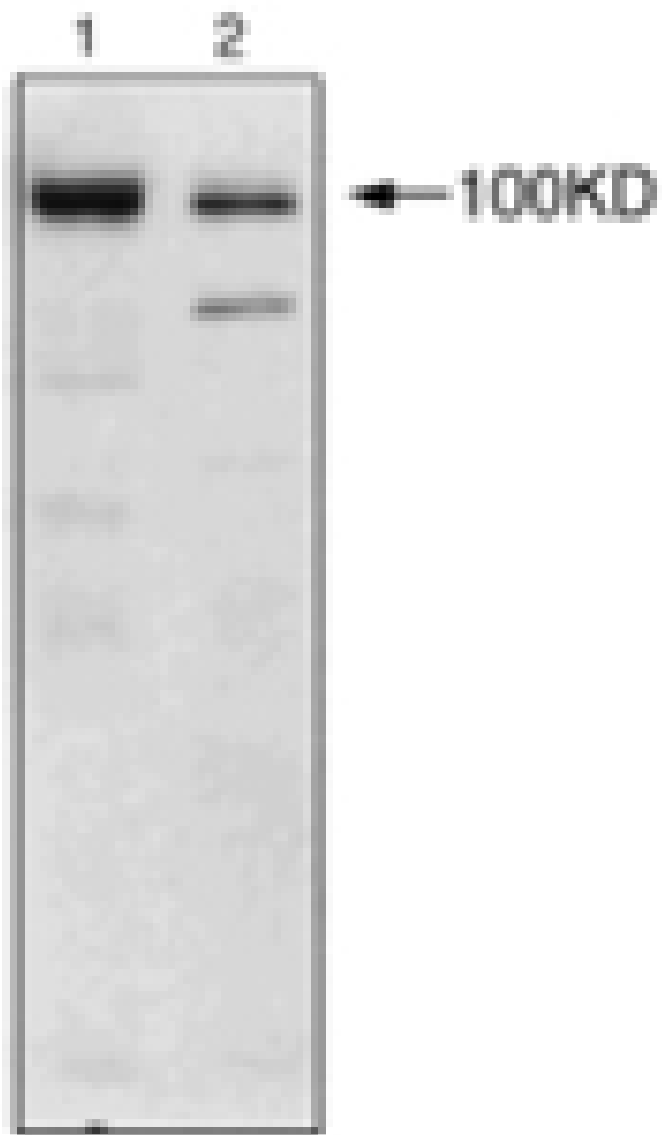
Ordering Information

[Order Online »](#)

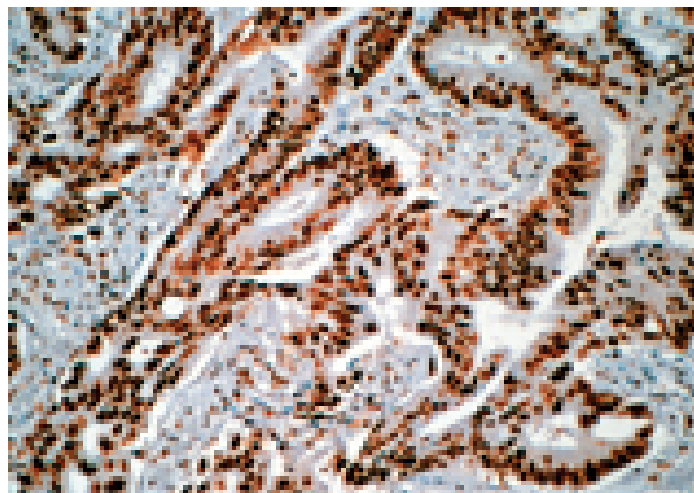
ADI-KAM-CP100-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HL60 (1) and ZR-75-1 (2) cell extracts, probed with Nucleolin mAb (4E2).



Immunohistochemistry analysis of nuclear staining of colon cancer sections using Nucleolin mAb (4E2).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Flow Cytometry, ICC, IHC, WB

Application Notes Detects a band of ~100kDa by Western blot.

Clone 4E2

Formulation Liquid. In PBS containing 50% glycerol.

Host Mouse

Immunogen Human nucleolin from Raji cell extract.

Isotype IgG1

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (5-10µg/ml final concentration)Immunocytochemistry (1-10µg/ml)Immunohistochemistry (1-10µg/ml)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P19338

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.948.9120
info@enzolifesciences.com
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com