

Nuclear pore complex proteins monoclonal antibody (MAb414) (purified)

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

Ordering Information

[Order Online »](#)

ENZ-ABS105-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

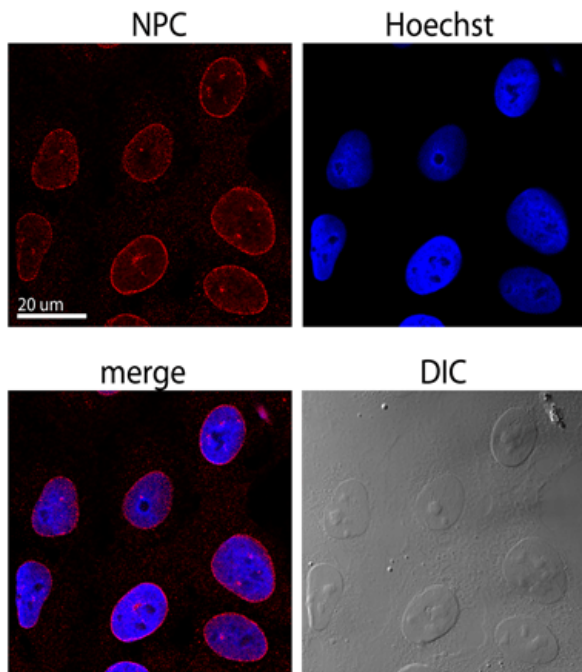


Figure: Human osteosarcoma cells stained with Nuclear pore complex proteins, mAb (MAb414) (purified) (Prod.no. ENZ-ABS105). Dilution 1:1000. Clockwise from top right you see NPC, Hoechst, merge, and differential interference contrast microscopy.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	NPC
Application	ICC, IHC, Immunoelectron microscopy, IP, WB
Clone	MAb414
Formulation	Liquid. In PBS containing 0.03% thimerosal.
Host	Mouse
Immunogen	Nuclear pore complex mixture.
Isotype	IgG1
Purity Detail	Protein G purified.
Species Reactivity	Vertebrate, Xenopus, Yeast
Specificity	Believed to recognize the repeated FXFG repeat sequence in nucleoporins including p62, p152 and p90 and other proteins.
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](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)