

Nuclear pore complex protein Nup153 monoclonal antibody (QE5)

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

Ordering Information

[Order Online »](#)

ENZ-ABS111-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Nucleoporin NUP153, NPC NUP153
Application	IF, IP, WB
Clone	QE5
Formulation	Liquid. In PBS containing 0.03% thimerosal.
Host	Mouse
Immunogen	Rat liver nuclear envelope proteins.
Isotype	IgG1
Purity Detail	Purified.
Species Reactivity	Eukaryote
Specificity	Recognizes eukaryote NUP153 as well as two related nuclear pore complex proteins: NUP214 and p62. By immunofluorescence, QE5 labels the nuclear envelope of eukaryotic cells giving a punctate staining pattern.
UniProt ID	P49791
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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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