

NeuN Recombinant monoclonal antibody (RM312)

RNA binding protein fox-1 homolog 3 (RBFOX3), also called Neuronal nuclei antigen and Fox-1 homolog C, is a pre-mRNA alternative splicing regulator. Regulates alternative splicing of RBFOX2 to enhance the production of mRNA species that are targeted for nonsense-mediated decay (NMD).

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

Ordering Information

[Order Online »](#)

ENZ-ABS671-0100	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	Neuronal nuclei antigen, RNA binding protein fox-1 homolog 3, Fox-1 homolog C, Fox-3
Application	IHC (PS), WB
Clone	RM312
Formulation	Liquid. In 1x PBS containing 50% glycerol, 1% BSA and 0.09% sodium azide
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the N-terminus of human NeuN/RBFOX3 .
Isotype	IgG
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (FFPE): 1:500-1:1,000Western Blot: 1:1,000-1:2,000Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from animal origin-free supernatant.
Species Reactivity	Human
UniProt ID	A6NFN3
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)