Neurofilaments (phosphorylated) monoclonal antibody (SMI31)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS215-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NF-H

Application ELISA, ICC, IHC, WB

Clone SMI-31

Formulation Liquid. In PBS containing thimerosal.

Host Mouse

lsotype IgG1

Purity Detail Purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:1,000)Western Blot

(1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Chicken, Human, Mammalian, Xenopus

Specificity Recognizes a phosphorylated epitope in extensively

phosphorylated neurofilament H and, to a lesser extent, in

neurofilament M.

UniProt ID P12036 (human)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

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

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