

NSF monoclonal antibody (9G7-3)

N-ethylmaleimide-sensitive factor (NSF) is involved in a variety of membrane fusion events and plays a prominent role in synaptic vesicle fusion. In order to interact with target membranes, NSF requires another set of soluble proteins called Soluble NSF attachment proteins (SNAPs). Together with synaptotagmin and alpha-SNAP, NSF modulates the interaction of SNAP25, VAMP, and syntaxin in an ATP-dependent manner to form a 20S complex involved in synaptic vesicle docking and fusion.

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

Citations: 2

[View Online »](#)

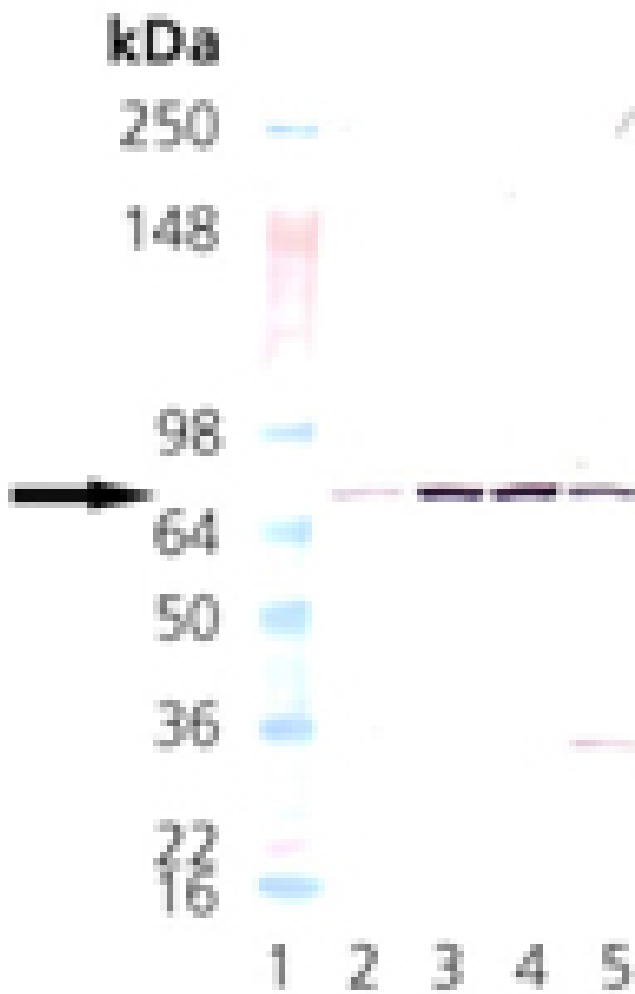
Ordering Information

[Order Online »](#)

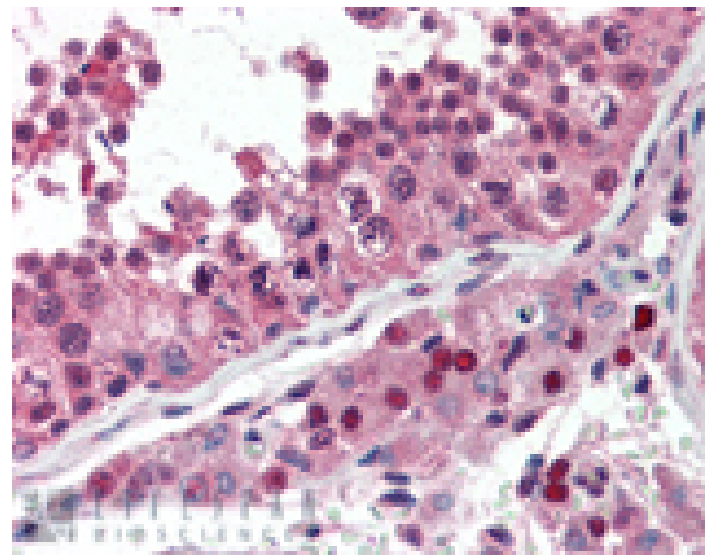
ADI-VAM-SV020-D	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: MW marker, Lane 2: HeLa (Heat Shocked), Lane 3: Vero, Lane 4: CHO-K1, Lane 5: ESK-4.



Immunohistochemistry analysis of human testis tissue stained with NSF, mAb (9G7-3) at 20 µg/ml.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-Ethylmaleimide-sensitive fusion protein

Application IHC (PS), WB

Application Notes Detects a band of ~76kDa by Western blot.

Clone 9G7-3

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Host Mouse

Immunogen Recombinant bovine NSF.

Isotype IgG2b

Purity Detail Protein G affinity purified.

Recommendation Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available
Dilutions/Conditions for all applications. Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Bovine, Chicken, Dog, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep

UniProt ID P46459 (human)



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 1 46 42 44 65
info-fr@enzolifesciences.com

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

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

Worry Free Guarantee This antibody is covered by our **Worry Free Guarantee**