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

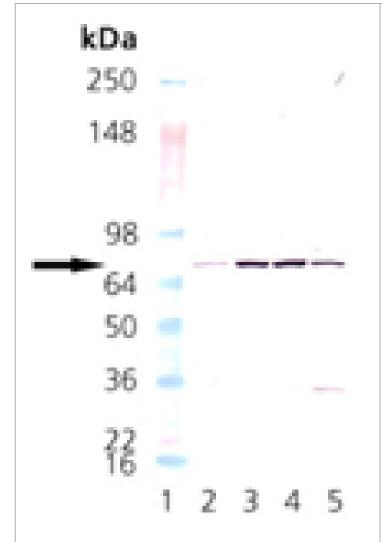
Order Online »

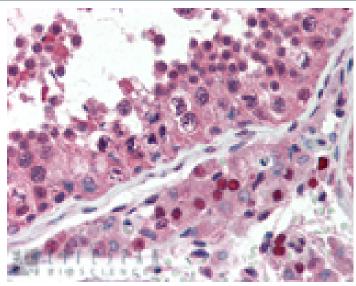
ADI-VAM-SV020-D

50µg

Manuals, SDS & CofA

View Online »





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

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

Lane 5: ESK-4.

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.

9G7-3 Clone

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

sodium azide.

Host Mouse

Immunogen Recombinant bovine NSF.

Isotype lgG2b

Purity Detail Protein G affinity purified.

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

> dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Purified from ascites. Source

Species Reactivity Bovine, Chicken, Dog, Hamster, Human, Monkey, Mouse,

Porcine, Rabbit, Rat, Sheep

UniProt ID P46459 (human)

Last modified: May 29, 2024



info-

eu@enzolifesciences.com