

Rabies Monoclonal Mouse Antibody (Rab-50)



Product Description

Reacts with viral glycoprotein of rabies virus strains SAD-Vnukovo and Pitman-Moore. This MAb is useful in detecting rabies virus by ELISA and Western. It is capable of neutralizing rabies virus. Rabies virus (Neurotropic virus) is a member of the Rhabdoviridae family. Rabies is a single stranded, neurotropic, negative sense RNA virus which encodes 5 proteins: a glycoprotein, a nucleoprotein, and three others. The mature virus has a bullet shape, a protein coat, and a lipid envelope. The outer surface of the virus is covered with thumb like glycoprotein projections 5-10 nm long and 3 nm in diameter. The virus averages approximately 780 nm in length. Lipid solvents destroy virus infectivity. Rabies virus is a very successful virus, with a very wide range of hosts. It causes an acute, central nervous system infection, characterized by CNS irritation, followed by paralysis and death. Approximately 50,000 human deaths each year are caused by rabies. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0943, Anti-Rabies (Rab-50)**

Product attributes	
Antibody number	#0943
Antibody reactivity (target)	Rabies
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	Rab-50
Isotype	IgG2b, kappa
Molecular weight	Not Known
Synonyms	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Purified virions of SAD-Vnukovo strain followed by Pitman-Moore strain.
Species reactivity	Rabies
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Rabies virus
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Endocrinology

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
BNCB	Biotin	N/A	N/A	N/A
BNUB	Purified	N/A	N/A	N/A
BNUM	Purified, BSA-free	N/A	N/A	N/A

Dye Features
[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.