Neurofilament H+L Monoclonal Mouse Antibody (2F11)



Product Description

This MAb reacts with a 200 kDa and 68 kDa protein, identified as heavy and light sub-units of neurofilaments (NF-H NF-L). Neurofilaments make up the main structural elements of axons and dendrites and are found in neurons, peripheral nerves, and sympathetic ganglion cells. Neurofilaments consist of three major subunits with molecular weights of 68 kDa (NF-L), 160 kDa (NF-M) and 200 kDa (NF-H). Anti-neurofilament stains a number of neural, neuroendocrine, and endocrine tumors. Neuromas, ganglioneuromas, gangliogliomas, ganglioneuroblastomas, and neuroblastomas stain positively for anti-neurofilament. Neurofilaments are also present in paragangliomas as well as adrenal and extra-adrenal pheochromocytomas. Carcinoids, neuroendocrine carcinomas of the skin, and oat cell carcinomas of the lung also express neurofilament. 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.

Call us: 800-304-5357

Product attributes				
Antibody number	#3912			
Antibody reactivity	Neurofilament H+L			
(target) Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	2F11			
Isotype	IgG1, kappa			
Molecular weight	200 kDa & 68 kDa			
Synonyms	NEFH; Neurofilament H; Neurofilament Heavy Polypeptide 200kDa; Neurofilament Triplet H Protein; NF-H; NF200. 68kDa neurofilament protein; Light molecular weight neurofilament protein; NEFL; Neurofilament light polypeptide 68kDa; Neurofilament light polypeptide; Neurofilament protein; light chain; Neurofilament subunit NF-L; Neurofilament triplet L protein; NF-L; NF-68			
Human gene symbol	NEFH & NEFL			
Entrez gene ID	4744			
SwissProt	P12036			
Unigene	198760			
Immunogen	Human NF-H from isolated brain cells			
Verified antibody	IHC (FFPE) (verified), WB (verified)			
Antibody target cellular localization Species reactivity	Cytoskeleton			
	Human, Mouse, Rat			
Positive control	IMR-32 cells. Cerebellum or Neuroblastoma.			
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			
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. mg/mL in PBS/0.05% BSA, Purified: 0. mg/mL in PBS/0.05% BSA/0.05% azid Purified, BSA-free: 1 mg/mL in PBS without azide			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
Cell/tissue expression	Neuroendocrine cells, Neurons			
Antibody research areas	Cancer, Cytoskeleton, Neuroscience			
Tumor expression	Lung cancer, Neuroendocrine cancer			

Email: btinfo@biotium.com

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

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.