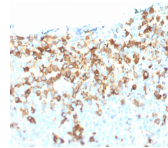


NGF Receptor Monoclonal Mouse Antibody (NGFR/1964)



Product Description

This antibody recognizes a glycoprotein of 75 kDa, identified as low affinity Nerve Growth Factor (NGF) Receptor (p75NGFR) or Neurotrophin Receptor (p75NTR). NGFR is expressed in various neural crest cells and their tumors such as melanocytes, melanomas, neuroblastomas, pheochromocytomas and neurofibromas. Reportedly, anti-NGFR is a reliable marker for desmoplastic and neurotropic melanomas. NGFR is expressed in mature non-neural cells such as perivascular cells, dental pulp cells, lymphoid follicular dendritic cells, basal epithelium of oral mucosa and hair follicles, prostate basal cells, and myoepithelial cells. Anti-NGFR stains the myoepithelial cells of breast ducts and intra-lobular fibroblasts of breast ducts.

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.

Product attributes

Antibody number	#1964
Antibody reactivity (target)	CD271, NGF-Receptor
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	NGFR/1964
Isotype	IgG2b, kappa
Molecular weight	75 kDa
Synonyms	CD271; Gp80-LNGFR; Low affinity nerve growth factor receptor; Low affinity neurotrophin receptor p75NTR; Nerve growth factor receptor (NGFR); p75 ICD; p75 Neurotrophin receptor; Tumor necrosis factor receptor superfamily member 16 (TNFRSF16)
Human gene symbol	NGFR
Entrez gene ID	4804
SwissProt	P08138
Unigene	415768 & 681726
Immunogen	Recombinant fragment of human p75 NGFR protein (around aa 281-421) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
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
Validated in protein array	Monospecific
Antibody research areas	Cancer, Developmental biology, Signal transduction
Cell/tissue expression	Neural crest cells
Tumor expression	Melanoma, Neuroendocrine cancer

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, BSA-free	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, of LI-COR Bioscience.

This datasheet was generated on August 31, 2024 at 12:15:52 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ngf-receptor-monoclonal-mouse-antibody-ngfr-1964/>