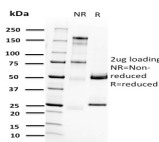


RET Proto-oncogene Monoclonal Mouse Antibody (RET/2662)



Product Description

The Ret proto-oncogene is structurally related to the growing family of tyrosine kinase transmembrane receptors and is involved in GDNF signaling. RET expression is reported in several regions of the central nervous system; in the developing cranial nerve ganglia and a subset of cells within dorsal root ganglia, in motor neurons in the spinal cord and hindbrain, in neuro-retina and the growing tips of the renal collecting ducts in developing kidney. Alterations in RET gene are associated with diseases including papillary thyroid carcinoma, multiple endocrine neoplasia (type 2A and 2B), familial medullary thyroid carcinoma, and a congenital developmental disorder known as Hirschsprung's disease.

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 2662, Anti-RET Proto-oncogene (RET/2662)

Product attributes

Antibody number	#2662
Antibody reactivity (target)	RET Proto-oncogene
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	RET/2662
Isotype	IgG2b, kappa
Molecular weight	150 kDa (precursor); 170 kDa (Mature)
Synonyms	Cadherin family member 12 (CDHF12); Cadherin related family member 16 (CDHR16); ELKS; HSCR1; Hydroxyaryl protein kinase; MEN2A; MEN2B; MTC1; Multiple endocrine neoplasia and medullary thyroid carcinoma 1; Oncogene RET; Proto-oncogene c-Ret; Proto-oncogene tyrosine-protein kinase receptor ret; PTC; RET; RET ELE1; RET51; RET9
Human gene symbol	RET
Entrez gene ID	5979
SwissProt	P07949
Unigene	350321
Immunogen	Recombinant fragment (around aa 702-848) of human RET protein (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane, Nucleus
Species reactivity	Human
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	Breast, Prostate or Colon Carcinoma.
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: 0.1 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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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
BNC74	CF®740	742/767	633-685	775/50	CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 26, 2024 at 04:22:43 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ret-proto-oncogene-monoclonal-mouse-antibody-ret-2662/>