

Renalase Monoclonal Mouse Antibody (RNLS/1940)



Product Description

Renalase is a 342 amino acid FAD-dependent amine oxidase that is highly expressed in kidney and is expressed at a lower level in heart, skeletal muscle and small intestine. Renalase is secreted in the blood by the kidney and it is thought to regulate cardiac function and systemic blood pressure. It is also suggested that Renalase functions as a hormone that metabolizes circulating catecholamines, which have an active role in the sympathetic and parasympathetic nervous systems. Individuals with chronic kidney disease and end-stage renal disease have markedly reduced levels of plasma Renalase than healthy individuals. Infusion of Renalase in animal models causes decrease in heart rate, cardiac contractility and blood pressure. Two isoforms of Renalase exist due to alternative splicing events.

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 1940, Anti-Renalase (RNLS/1940)

Product attributes

Antibody number	#1940
Antibody reactivity (target)	Renalase
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	RNLS/1940
Isotype	IgG2b, kappa
Molecular weight of antigen	38 kDa
Synonyms	Chromosome 10 open reading frame 59; MAO C; MAO-C; Monoamine oxidase-C; Renalase; Renalase FAD dependent amine oxidase; RNLS
Human gene symbol	RNLS
Entrez gene ID	55328
SwissProt	Q5VYX0
Unigene	149849
Immunogen	Recombinant human RNLS protein fragment (around aa 34-235) (exact sequence is proprietary)
Antibody target cellular localization	Secreted (extracellular)
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	293T Cells. Heart or Kidney
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, Metabolism
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 June 13, 2026 at 08:00:31 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/renalase-monoclonal-mouse-antibody-rnls-1940/>