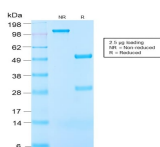


IGF-1 Recombinant Monoclonal Rabbit Antibody (IGF1/2872R)



Product Description

This antibody is specific to Insulin-like Growth Factor (IGF-1) and shows minimal cross-reaction with IGF-11, Proinsulin, MSF, and Insulin. IGF-1 is a polypeptide growth factor with two isoforms that are produced by alternative splicing. Isoform 1 is also known as IGF-1B while isoform 2 is known as IGF-1A. IGF-1 stimulates the proliferation of a wide range of cell types including muscle, bone and cartilage tissue. It functions as an autocrine regulator of growth. Activation of IGF system has emerged as a key factor for tumor progression and resistance to apoptosis in many cancers like those of breast, thyroid and colon.

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 2872, Anti-Insulin-like Growth Factor-1 (IGF1/2872R)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2872
Antibody reactivity (target)	Insulin-like Growth Factor-1
Antibody type	Primary
Host species	Rabbit
Clonality	Monoclonal, Recombinant Monoclonal
Clone	IGF1/2872R
Isotype	IgG
Molecular weight of antigen	~7.6 kDa
Synonyms	Insulin-like Growth Factor IA (IGF-1A); Insulin-like Growth Factor IB (IGF-1B); Mechano Growth Factor (MGF); Somatomedin-C
Human gene symbol	IGF1
Entrez gene ID	3479
SwissProt	P05019
Unigene	160562
Immunogen	Recombinant full-length human IGF-1 protein
Antibody target cellular localization	Secreted (extracellular)
Verified antibody applications	Flow (verified)
Species reactivity	Human, Mouse, Rabbit, Rat
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL, ELISA For coating, order Ab without BSA, Optimal dilution for a specific application should be determined by user
Positive control	Pancreas or brain. Breast, Thyroid or Colon Cancers; IGF-1 recombinant protein
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	Cancer, Metabolism
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Tumor expression	Breast cancer, Colorectal cancer, Thyroid 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
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 8, 2026 at 02:37:39 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/insulin-like-growth-factor-1-recombinant-monoclonal-rabbit-antibody-igf1-2872r/>