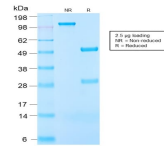


# IGF-1 Recombinant Monoclonal Rabbit Antibody (IGF1/2872R)



## 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-IB while isoform 2 is known as IGF-IA. 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

## Product attributes

<b>Antibody number</b>	#2872
<b>Antibody reactivity (target)</b>	Insulin-like Growth Factor-1
<b>Antibody type</b>	Primary
<b>Host species</b>	Rabbit
<b>Clonality</b>	Recombinant Monoclonal
<b>Clone</b>	IGF1/2872R
<b>Isotype</b>	IgG
<b>Molecular weight</b>	~7.6 kDa
<b>Synonyms</b>	Insulin-like Growth Factor IA (IGF-IA); Insulin-like Growth Factor IB (IGF-IB); Mechano Growth Factor (MGF); Somatomedin-C
<b>Human gene symbol</b>	IGF1
<b>Entrez gene ID</b>	3479
<b>SwissProt</b>	P05019
<b>Unigene</b>	160562
<b>Immunogen</b>	Recombinant full-length human IGF-1 protein
<b>Antibody target cellular</b>	Secreted (extracellular)
<b>Verified antibody applications</b>	Flow (verified)
<b>Species reactivity</b>	Human, Mouse, Rabbit, Rat
<b>Antibody application notes</b>	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
<b>Positive control</b>	Pancreas or brain, Breast, Thyroid or Colon Cancers; IGF-1 recombinant protein
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer, Metabolism
<b>Tumor expression</b>	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	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
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.