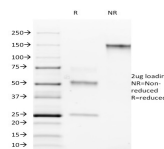


# ROR-Gamma / RORC Monoclonal Mouse Antibody (RORC/2942)



## Product Description

This MAbs recognizes a protein of 63 kDa, identified as ROR-C. Its epitope maps in between aa1-50. The nuclear orphan receptors ROR and ROR are members of the nuclear hormone receptor superfamily. Members of this family act by directly associating with DNA sequences known as hormone response elements (HREs) and typically bind DNA as either homo- or heterodimers. RORalpha and RORgamma are unique in that they bind DNA as monomers. RORalpha has multiple isoforms that share common DNA and putative ligand-binding domains, but differ in their amino terminal domains, which are generated by alternative RNA processing. RORgamma comprises a 560 amino acid protein that shares 50% amino acid identity with RORalpha and is most highly expressed in skeletal muscle. Although these proteins are considered orphan receptors, due to a lack of defined ligands, experimental evidence has shown that melatonin may be the natural ligand for these nuclear receptors.

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.

**Catalog number key for antibody number 2942, Anti-ROR-Gamma|RORC (RORC/2942)**

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

## Product attributes

Antibody number	#2942
Antibody reactivity (target)	ROR-Gamma, RORC
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	RORC/2942
Isotype	IgG2a, kappa
Molecular weight	63 kDa
Synonyms	Nuclear receptor ROR-gamma; Nuclear receptor RZR-gamma; Nuclear receptor subfamily 1 group F member 3 (NR1F3); RAR related orphan receptor C; Retinoid-related orphan receptor-gamma (RORG); Rorc; RZR GAMMA; RZRG; TOR
Human gene symbol	RORC
Entrez gene ID	6097
SwissProt	P51449
Unigene	256022; 607993
Immunogen	Recombinant full-length human RORC protein
Antibody target cellular localization	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	Human Lymphocytes. Liver or skeletal muscle tissue.
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 (Bos taurus), 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	<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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on December 13, 2025 at 08:36:01 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/ror-gamma-rorc-monoclonal-mouse-antibody-rorc-2942/>