FCGRT / FcRn Monoclonal Mouse Antibody (FCGRT/2932)



Product Description

This gene encodes a receptor that binds the Fc region of monomeric immunoglobulin G. FCGRT is present in the intestinal epithelium of neonates and mediates the selective uptake of immunoglobulin G (IgG) in mothers' milk, thereby helping newborn to acquire passive immunity. FCGRT is comprised of a heavy chain and -2-Microglobulin. FCGRT localizes in endosomes of vascular endothelial cells and selectively recycles IgG to the cell surface, thus protecting IgG from lysosomal degradation.

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 2932, Anti-FCGRT/FcRn (FCGRT/2932)

Call us: 800-304-5357 Email: techsupport@biotium.com

Product attributes			
Antibody number	#2932		
Antibody reactivity (target)	FCGRT, FcRn		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	FCGRT/2932		
Isotype	IgG2b, kappa		
Molecular weight	46 kDa		
Synonyms	Fc fragment of IgG; receptor transporter; alpha; FCGRT; FCRN alpha chain; IgG Fc fragment receptor transporter alpha chain; IgG Gc receptor; IgG receptor FcRn large subunit p51; Immunoglobulin receptor; intestinal; heavy chain; Neonatal Fc receptor		
Human gene symbol	FCGRT		
Entrez gene ID	2217		
SwissProt	P55899		
Unigene	111903		
Immunogen	Recombinant fragment of human FCGRT protein (around aa 24-215) (Exact sequence is proprietary)		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Positive control	THP1 or HepG2 cells. Cerebellum or Testis.		
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/in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/m 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 Chipage hamster every calls lamite for the prodific let		

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 August 20, 2025 at 10:48:59 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fcgrt-fcrn-monoclonal-mouse-antibody-fcgrt-2932/