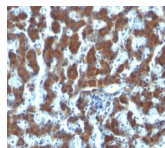


Retinol Binding Protein 1 Monoclonal Mouse Antibody (G4E4)



Product Description

This MAb recognizes an epitope within the 74-182 C-terminal sequence (11kD peptide fragment) of human serum Cellular Retinol Binding Protein 1 (CRBP 1), a single-chain glycoprotein belonging to the superfamily of hydrophobic molecule transporter proteins, which is responsible for transport of retinol (vitamin A1) from the liver to peripheral target tissues, like the eye, where it mediates the cellular uptake. CRBP 1 is synthesized by hepatic parenchymal cells where it becomes bound to its ligand retinol and is then released into the circulation, where it binds further to the protein transthyretin, to form a transporting complex, which is big enough not to be lost by filtration through the kidney glomeruli. It is detected in nearly all tissues with higher expression in adult ovary, pancreas, pituitary gland, adrenal gland, and fetal liver.

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 0855, Anti-Retinol Binding Protein 1 (G4E4)

Product attributes

Antibody number	#0855
Antibody reactivity (target)	Retinol Binding Protein 1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	G4E4
Isotype	IgG1, kappa
Molecular weight	21-25 kDa
Synonyms	Cellular retinol-binding protein 1, CRBP1, CRBP2, RBP1, RBP2, RBP4, RBPC, Retinol binding protein 1, Retinol binding protein 1 cellular, Retinol binding protein 2 cellular, Retinol binding protein 4 plasma
Human gene symbol	RBP1
Entrez gene ID	5947, 5948 & 5950
SwissProt	P02753, P09455 & P50120
Unigene	529571
Immunogen	Retinol binding protein-1 purified from human plasma
Antibody target cellular location	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Chimpanzee, Goat, Human, Monkey, 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. Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user
Positive control	Hepatic or ovarian carcinoma
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	Metabolism

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)

[CF®488A Features](#)

[CF®568 Features](#)

[CF®594 Features](#)

[CF®640R Features](#)

[CF®647 Features](#)

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, or of LI-COR Bioscience.