## Retinol Binding Protein 1 Monoclonal Mouse Antibody (RBP/872)



## **Product Description**

Recognizes a protein of 21 kDa-25 kDa, identified as retinol binding protein-1 (RBP1). This protein belongs to the lipocalin family and is the specific carrier for retinol (vitamin A alcohol) in the blood. It delivers retinol from the liver stores to the peripheral tissues. In plasma, the RBP-retinol complex interacts with transthyretin, which prevents its loss by filtration through the kidney glomeruli. A deficiency of vitamin A blocks secretion of the binding protein post-transnationally and results in defective delivery and supply to the epidermal cells.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0872, Anti-Retinol Binding Protein 1 (RBP/872)

## Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#0872			
Antibody reactivity (target)	Retinol Binding Protein 1			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	RBP/872			
Isotype	IgG1, kappa			
Molecular weight	21-25 kDa			
Synonyms	Cellular retinol-binding protein I, 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	Recombinant human retinol binding protein-1 (RBP1)			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Secreted (extracellular)			
Species reactivity	Goat, Human, Monkey, Mouse, Rabbit, Rat			
Antibody application notes	Immunohistology formalin-fixed 1-2 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, immunofluorescence 1-2 ug/mL, 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 i PBS without azide			
Antibody research areas	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.			

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 November 22, 2025 at 11:08:43 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-retinol-binding-protein-1-rbp1872/">https://biotium.com/product/monoclonal-anti-retinol-binding-protein-1-rbp1872/</a>