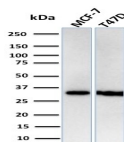


Replication Protein A2 Recombinant Monoclonal Rabbit Antibody (RPA2/3140R)



Product Description

Recognizes a protein of 34 kDa, identified as a subunit of Replication Protein A (RPA) (also known as human single-stranded DNA binding protein, or HSSB). RPA from human cells is a stable heterotrimer of 70 kDa, 32-34 kDa, and 11-14 kDa subunits (RPA70, RPA32, and RPA14 respectively). It is involved in DNA replication, repair, and recombination. RPA is required for the SV40 large tumor antigen-catalyzed unwinding of SV40 DNA and stimulates DNA polymerase (pol) alpha and delta. RPA34 is phosphorylated at the G1/S boundary of the cell cycle or upon exposure of cells to DNA damage-inducing agents including ionizing and UV radiation.

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 3140, Anti-Replication Protein A2 (RPA2/3140R)

Product attributes

Antibody number	#3140
Antibody reactivity (target)	Replication Protein A2
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	RPA2/3140R
Isotype	IgG
Molecular weight	32-34 kDa
Synonyms	p32; p34; REPA2; Replication protein A 32kDa subunit; Replication protein A 34kDa subunit; Replication protein A2; RF-A protein 2; Rf-A2; RFA; RP-A p32; RP-A p34; RP21C; RPA32; RPA34; RPLP1; RPLP2
Human gene symbol	RPA2
Entrez gene ID	6118
SwissProt	P15927
Unigene	703070; 79411
Immunogen	Recombinant human full-length Replication protein A2 (RPA2) protein
Verified antibody applications	WB (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined.
Positive control	MCF-7, HeLa or HepG2 cell lysates (WB). Human tonsil (IHC).
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
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	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 December 12, 2025 at 02:15:12 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/replication-protein-a2-recombinant-monoclonal-rabbit-antibody-rpa2-3140r/>