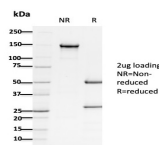


Adenomatous Polyposis Coli / APC Monoclonal Mouse Antibody (ALi 12-28)



Product Description

Adenomatous Polyposis Coli (APC) is a protein with many cellular functions. It is a tumor suppressor and negative regulator of Wnt signaling. APC has roles in cell adhesion and cell migration. Mutations in APC are associated with several diseases. The adenomatous polyposis syndromes, familial adenomatous polyposis (FAP) and Gardner's syndrome (GS), are characterized by numerous adenomatous polyps throughout the entire colon. These polyps invariably progress to colon cancer in addition to other extracolonic manifestations.

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 3076, APC / Adenomatous Polyposis Coli / FAP Monoclonal Mouse Antibody (ALi 12-28)

Product attributes

research-areas	Cytoskeleton, Signal transduction
Antibody number	#3076
Antibody reactivity (target)	Adenomatous Polyposis Coli
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ALi 12-28
Isotype	IgG1, kappa
Molecular weight	66 kDa
Synonyms	Adenomatous Polyposis Coli; APC; CC1; Deleted in polyposis 2.5; Protein Phosphatase 1, Regulatory Subunit 46; DP2; DP3
Human gene symbol	APC
Entrez gene ID	324
SwissProt	P25054
Unigene	158932
Immunogen	Recombinant fragment of human APC protein (around aa 1-433)
Antibody target cellular localization	Golgi apparatus, Plasma membrane, Nucleus
Species reactivity	Human
Expected antibody applications	IHC (frozen) (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
Positive control	COLO 320DM whole cell lysate. HCT116 cells. Tonsil, placenta, breast or colon.
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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- PNAS USA (1998) 95:3122-3127. (WB, IF, epitope mapping)
- EMBO Rep (2005) 6:184-190. (IF)
- BMC Cell Biol (2006) 7: 3. (IF, WB)
- Int J Mol Med (2008) 21:19-31. (IHC, frozen)
- J Biol Chem (2012) 287(34): 28552-28563. (IP, WB)

This datasheet was generated on September 11, 2025 at 01:35:50 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/apc-adenomatous-polyposis-coli-fap-monoclonal-mouse-antibody-ali-12-28/>