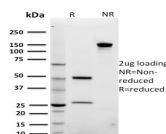


ADP-Ribosylation Factor 1 (ARF1) Monoclonal Mouse Antibody (3F1)



Product Description

The ADP-ribosylation factor (ARF) family comprises a group of structurally and functionally conserved proteins, which are members of the Ras superfamily of regulatory GTP-binding proteins. The ARF family is divided functionally into the ARF and the ARF-like proteins. ARF's share more than 60% sequence identity, appear to be ubiquitous in eukaryotes, and are highly conserved evolutionarily. ARF is involved in intracellular protein traffic to and within the Golgi complex. ARF has a number of disparate activities including maintenance of organelle integrity, assembly of coat proteins, as a co-factor for cholera toxin and as an activator of phospholipase D.

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 2029, ARF1 Monoclonal Mouse Antibody (3F1)

Product attributes

Antibody number	#2029
research-areas	Membrane trafficking, Organelle markers
Antibody reactivity (target)	ADP-Ribosylation Factor 1, ARF1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3F1
Isotype	IgG2a, kappa
Molecular weight	20 kDa
Synonyms	ADP-ribosylation factor 1; ARF1
Human gene symbol	ARF1
Entrez gene ID	375
SwissProt	P84077
Unigene	286221
Immunogen	Synthetic peptide SNQLRNQ, corresponding to C terminal amino acids 174-180 of Human ARF1.
Antibody target cellular localization	Golgi apparatus
Species reactivity	Human, Mouse, Rat
Expected antibody applications	IF (published for clone), WB (published for clone)
Positive control	HeLa cells. Kidney, Liver, Brain or Breast.
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.

1. J Biol Chem (2009) 284(20): 13948-13957. (WB)
2. J Biochem (2016) 159(2): 201-208. (IF; WB)

This datasheet was generated on August 4, 2025 at 04:50:33 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/arf1-monoclonal-mouse-antibody-3f1/>