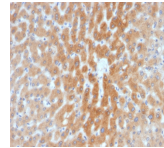


# ADP-Ribosylation Factor 1 (ARF1) Monoclonal Mouse Antibody (1A9/5)



## 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 2028, ARF1 Monoclonal Mouse Antibody (1A9/5)**

## Product attributes

Antibody number	#2028
research-areas	Membrane trafficking, Organelle markers
Antibody reactivity (species)	ADP-Ribosylation Factor 1, ARF1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	1A9/5
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	C-terminal human ARF1 peptide SNQLRNQ
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Golgi apparatus
Species reactivity	Human, Mouse, Rat
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

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