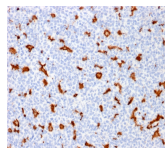


Allograft Inflammatory Factor 1 / IBA1 Monoclonal Mouse Antibody (AIF1/1909)



Product Description

Allograft Inflammatory Factor 1 (also known as AIF1 or IBA1) is a cytoplasmic, calcium-binding protein that is thought to play a role in macrophage activation and function. AIF1, containing two EF domains, is induced by cytokines and Interferons. In an unstimulated state, AIF1 colocalizes with actin, and upon stimulation, translocates to lamellipodia. It is also a marker of human microglia and is expressed by macrophages in injured skeletal muscle. The gene encoding AIF1 resides in the tumor necrosis factor (TNF) cluster of genes, located in the region represented by the human major histocompatibility complex (MHC). 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 1909, Anti-AIF1 / Iba1 (AIF1/1909)**

Product attributes

research-areas	Immunology
Antibody number	#1909
Antibody reactivity (species)	Allograft Inflammatory Factor 1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	AIF1/1909
Isotype	IgG1, kappa
Molecular weight	17 kDa
Synonyms	Allograft inflammatory factor 1; AIF1; IBA1
Human gene symbol	AIF1
Entrez gene ID	199
SwissProt	P55008
Unigene	76364
Immunogen	Purified fragment of human recombinant AIF1 protein (around aa 1-146) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	Lymph Node, Tonsil or Kidney
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 in PBS without azide
Validated in protein array	Monospecific
Antibody research areas	Immunology
Cell/tissue expression	Microglia, Monocytes/macrophages

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
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 LI-COR Bioscience.