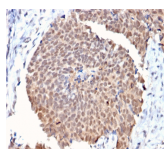


Fas associated factor 1 Monoclonal Mouse Antibody (CPTC-FAF1-2)



Product Description

In contrast to growth factors which promote cell proliferation, FAS ligand (FAS-L) and the tumor necrosis factors (TNFs) rapidly induce apoptosis. Cellular response to FAS-L and TNF is mediated by structurally related receptors containing a conserved 'death domain' and belonging to the TNF receptor superfamily. TRADD, FADD and RIP are FAS/TNF-RI interacting proteins that contain a death domain homologous region (DDH). TRADD (TNF-RI-associated death domain) and FADD (FAS-associated death domain) associate with the death domains of both FAS and TNF-RI via their DDH regions, while RIP associates exclusively with FAS. An additional FAS interacting protein designated FAF1, for FAS-associated protein factor-1, binds with the cytoplasmic tail of wildtype but not LPR mutant FAS. When overexpressed in cells, FAF1 enhances the efficiency of FAS-mediated apoptosis. In contrast to TRADD, FADD and RIP, FAF1 lacks a DDH and cannot induce apoptosis independently of FAS activation.

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 2268, Anti-Fas associated factor 1 (CPTC-FAF1-2)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2268
Antibody reactivity (target)	Fas associated factor 1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-FAF1-2
Isotype	IgG2b, kappa
Molecular weight	75-80 kDa
Synonyms	CGI-03; FAS-associated factor 1; HFAF1s; UBX domain-containing protein 12; UBX domain-containing protein 3A
Human gene symbol	FAF1
Entrez gene ID	11124
SwissProt	Q9UNN5
Unigene	530402
Immunogen	Recombinant human full-length FAF1 protein
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Positive control	HEK293, MCF-7, HeLa whole cell lysate. Human testis, prostate, breast carcinoma or colon carcinoma.
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
Validated in protein array	Monospecific
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 03:15:24 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/fas-associated-factor-1-monoclonal-mouse-antibody-cptc-faf1-2/>