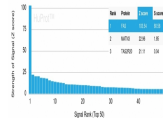


CD95 Monoclonal Mouse Antibody (FAS/3588)



Product Description

This MAbs specifically recognizes CD95, also known as Fas, a transmembrane glycoprotein with a MW of 40-45 kDa, containing 8 kDa of N-glycoside-linked polysaccharide. It is a receptor for TNFSF6/FASLG, a member of the nerve growth factor receptor/tumor necrosis factor superfamily, mediating receptor-triggered apoptosis. The adapter molecule FADD recruits caspase-8 to the activated receptor. The resulting death-inducing signaling complex (DISC) performs caspase-8 proteolytic activation, which initiates the subsequent cascade of caspases (aspartate-specific cysteine proteases) mediating apoptosis. FAS-mediated apoptosis may have a role in the induction of peripheral tolerance, in the antigen-stimulated suicide of mature T-cells, or both. The secreted isoforms 2 to 6 block apoptosis (in vitro). CD95 antigen is expressed on the surface of various cell types, preferentially on the CD45RA^{low} CD45RO^{high} subset of memory T lymphocytes.

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 3588, Anti-CD95|FAS|TNFRSF6 (FAS/3588)

Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

Product attributes

Antibody number	#3588
Antibody reactivity (target)	CD95, FAS, TNFRSF6
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FAS/3588
Isotype	IgG2b, kappa
Molecular weight	38-50 kDa
Synonyms	ALPS1A; Apoptosis APO1 antigen 1; Apoptosis-mediating surface antigen FAS; APT1; Delta Fas; Fas AMA; Fas cell surface death receptor; FAS1; FASL receptor; FASLG receptor; FASTM; Surface antigen APO1; Tumor necrosis factor receptor superfamily member 6 (TNFRSF6)
Human gene symbol	FAS
Entrez gene ID	355
SwissProt	P25445
Unigene	244139
Immunogen	Recombinant fragment (around aa26-96) of human CD95 (FAS) protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	MCF-7 cells (IF/FACS). Human hepatocellular or bladder carcinoma (IHC).
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 July 19, 2025 at 05:15:51 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd95-monoclonal-mouse-antibody-fas-3588/>