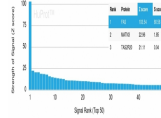


# CD95 Monoclonal Mouse Antibody (FAS/3588)



## Product Description

This MAb 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<sup>low</sup> CD45RO<sup>high</sup> subset of memory T lymphocytes. Primary antibodies are available purified, or with a selection of fluorescent CF<sup>®</sup> dyes and other labels. CF<sup>®</sup> dyes offer exceptional brightness and photostability. See the [CF<sup>®</sup> Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF<sup>®</sup>405S and CF<sup>®</sup>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<sup>®</sup> 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.

## Product attributes

<b>Antibody number</b>	#3588
<b>Antibody reactivity (target)</b>	CD95, FAS, TNFRSF6
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	FAS/3588
<b>Isotype</b>	IgG2b, kappa
<b>Molecular weight</b>	38-50 kDa
<b>Synonyms</b>	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)
<b>Human gene symbol</b>	FAS
<b>Entrez gene ID</b>	355
<b>SwissProt</b>	P25445
<b>Unigene</b>	244139
<b>Immunogen</b>	Recombinant fragment (around aa26-96) of human CD95 (FAS) protein (exact sequence is proprietary)
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Human
<b>Positive control</b>	MCF-7 cells (IF/FACS). Human hepatocellular or bladder carcinoma (IHC).
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Validated in protein assays</b>	Monospecific
<b>Shelf life</b>	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 <sup>®</sup>	<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.