Phospho C-Jun Monoclonal Mouse Antibody (C-J 4C4/1)



Product Description

This gene is the putative transforming gene of avian sarcoma virus 17. It encodes a protein which is highly similar to the viral protein, and which interacts directly with specific target DNA sequences to regulate gene expression. This gene is intronless and is mapped to 1p32-p31, a chromosomal region involved in both translocations and deletions in human malignancies.

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 2282, Anti-Phospho C-Jun (C-J 4C4/1)

Call us : <u>800-304-5357</u>

Product attributes			
Antibody number	#2282		
Antibody reactivity (target)	Phospho C-Jun		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	C-J 4C4/1		
Isotype	IgG1, kappa		
Molecular weight	39 kDa		
Synonyms	Activator protein 1 (API); cJun; Enhancer Binding Protein AP1; Jun Activation Domain Binding Protein; Jun proto oncogene; Transcription Factor AP1; V jun avian sarcoma virus 17 oncogene homolog; VV jun sarcoma virus 17 oncogene homolog; V-jun avian sarcoma virus 17 oncogene homolog; vJu Avian Sarcoma Virus 17 oncogene Homolog		
Human gene symbol	JUN		
Entrez gene ID	3725		
SwissProt	P05412		
Unigene	696684		
Immunogen	c-Jun protein phosphorylated at T91 and T93		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Nucleus		
Species reactivity	Human, Rat		
Positive control	HEK293, HeLA or A-431 cells; Human normal lung or bladder tissue.		
Shipping condition	Room temperature		
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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/r in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/ml PBS without azide		
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.		

Austinador # wrativ	Conjugation	Fx/Fxx (xxxx)	I agay lina	Detection channel	Due Feetures
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 November 21, 2025 at 06:28:00 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/phospho-c-jun-monoclonal-mouse-antibody-c-j-4c4-1/