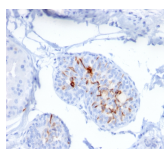


ZNF690 / ZSCAN29 Monoclonal Mouse Antibody (ZSCAN29/2610)



Product Description

Zinc finger proteins contain DNA-binding domains and have a wide variety of functions, most of which encompass some form of transcriptional activation or repression. The majority of zinc finger proteins contain a Kruppel-type DNA binding domain and a KRAB domain, which is thought to interact with KAP1, thereby recruiting histone modifying proteins. Zinc finger protein 690 (ZNF690), also known as ZSCAN29, is a 851 amino acid member of the Kruppel C2H2- type zinc finger protein family. Localized to the nucleus, ZNF690 contains six C2H2-type zinc fingers and one KRAB domain through which it is thought to be involved in DNA-binding and transcriptional regulation. Four isoforms of ZNF690 exist as a result of alternative splicing events.

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 2610, Anti-ZNF690/ZSCAN29 (ZSCAN29/2610)

Product attributes

Antibody number	#2610
Antibody reactivity (target)	ZNF690, ZSCAN29
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ZSCAN29/2610
Isotype	IgG1, kappa
Molecular weight of antigen	97 kDa
Synonyms	FLJ35867; KOX31 like zinc finger protein; MGC129894; MGC129895; zinc finger and SCAN domain-containing 29 (ZSCAN29); Zinc finger protein 690 (ZNF690)
Human gene symbol	ZSCAN29
Entrez gene ID	146050
SwissProt	Q8IYW8
Unigene	418287
Immunogen	Recombinant full-length human ZSCAN29 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human
Positive control	HeLa whole cell lysate. Tonsil. Breast 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 (<i>Bos taurus</i>), 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 June 4, 2026 at 02:46:17 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/znf690-zscan29-monoclonal-mouse-antibody-zscan29-2610/>