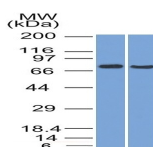


TCF4 Monoclonal Mouse Antibody (TCF4/1705)



Product Description

This antibody recognizes a protein of 71 kDa, identified as Transcription Factor 4 (TCF4). It is a basic helix-loop-helix transcription factor. The protein recognizes an Ephrussi-box (E-box) binding site (CANNTG) – a motif first identified in immunoglobulin enhancers. This gene is broadly expressed, and may play an important role in nervous system development. Defects in this gene are a cause of Pitt-Hopkins syndrome.

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 1705, Anti-TCF4 (TCF4/1705)

Product attributes

Antibody number	#1705
Antibody reactivity (target)	TCF4
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TCF4/1705
Isotype	IgG2a, kappa
Molecular weight	71 kDa
Synonyms	bHLHb19; Class B basic helix-loop-helix protein 19; Immunoglobulin transcription factor 2; ITF-2; PTHS; SEF2; SL3-3 enhancer factor 2; TCF4; Transcription factor 4
Human gene symbol	TCF4
Entrez gene ID	6925
SwissProt	P15884
Unigene	605153
Immunogen	Recombinant human TCF4 protein fragment (around aa 365-671) (exact sequence is proprietary)
Antibody target cellular localization	Nucleus
Verified antibody applications	WB (verified)
Species reactivity	Human
Antibody application notes	ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	HeLa or HepG2 cells. Bladder
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Developmental biology, Transcription factors
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 November 29, 2024 at 12:45:03 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/tcf4-monoclonal-mouse-antibody-tcf4-1705/>