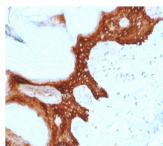


14-3-3 Sigma / Stratifin Monoclonal Mouse Antibody (CPTC-SFN-2)



Product Description

Stratifin is a p53-induced tumor suppressor gene which is activated in response to DNA damage, causing cell cycle arrest at G2 phase by blocking cdc2-cyclin B1 complex from entering the nucleus. It is inactivated in breast, lung, prostate, liver and gastric cancer. It is associated with poor prognosis when its down-regulation is observed in epithelial ovarian cancer. SFN expression could contribute to cancer cell proliferation and the development and/or progression of human gastrointestinal cancer.

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 2481, 14-3-3 Sigma / Stratifin Monoclonal Mouse Antibody (CPTC-SFN-2)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2481
Antibody reactivity (target)	14-3-3 Sigma, Stratifin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Synonyms	Stratifin; YWHAS
Clone	CPTC-SFN-2
Isotype	IgG1, kappa
Molecular weight	28 kDa
Human gene symbol	SFN
Entrez gene ID	2810
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
SwissProt	P31947
Unigene	523718
Immunogen	Recombinant human full-length SFN protein
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	PC-3 cells, Skin and tissues enriched in stratified squamous keratinizing epithelium. HaCaT, hTCEpi, A-431 cell lines.
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
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
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.
Regulatory status	For research use only (RUO)
Validated in protein array	Monospecific

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 December 27, 2025 at 09:46:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/14-3-3-sigma-stratifin-monoclonal-mouse-antibody-cptc-sfn-2/>