## TTF-1 / NKX2.1 Recombinant Monoclonal Rabbit Antibody (NX2.1/1855R)



## **Product Description**

Antibody # prefix

**BNUM** 

This antibody recognizes a protein of 40 kDa, identified as Thyroid transcription factor-1 (TTF-1). TTF-1 is a member of the NKx2 family of homeodomain transcription factors. It is expressed in epithelial cells of the thyroid gland and the lung. Nuclei from liver, stomach, pancreas, small intestine, colon, kidney, breast, skin, testes, pituitary, prostate, and adrenal glands are unreactive. Anti-TTF-1 is useful in differentiating primary adenocarcinoma of the lung from metastatic carcinomas originating in the breast, mediastinal germ cell tumors, and malignant mesothelioma. It can also be used to differentiate small cell lung carcinoma from lymphoid infiltrates. Loss of TTF-1 expression in non-small cell lung carcinoma has been associated with aggressive behavior of such neoplasms. TTF-1 reactivity is also seen in thyroid 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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1855, Anti-TTF-1 / NKX2.1 (NX2.1/1855R)

Conjugation

Purified,

BSA-free

## **Product attributes**

Fiduuci attiibutes				
Antibody number	#1855			
Antibody reactivity (target)	NKX2.1, TTF-1			
Antibody type	Primary			
Host species	Rabbit			
Clonality	Recombinant Monoclonal			
Clone	NX2.1/1855R			
Isotype	IgG			
Molecular weight	40 kDa			
Synonyms	BCH; Benign chorea; BHC; Homeobox protein Nkx2.1; NK2 homeobox 1; NKX2.1; NKX2A; TEBP; Thyroid nuclear factor-1; Thyroid specific enhancer binding protein; Thyroid transcription factor-1 (TTF-1); Tin man; TITF1; TTF-1			
Human gene symbol	NKX2-1			
Entrez gene ID	7080			
SwissProt	P43699			
Unigene	94367			
Immunogen	Recombinant human full-length TTF-1 protein			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Nucleus			
Species reactivity	Human, Mouse, Rat			
Positive control	MAD109, MLE-15, H441-4, or H345 cells. Normal thyroid or lung.			
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			
Antibody research areas	Cancer, Transcription factors			
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.			
Tumor expression	Lung cancer, Thyroid cancer			

Dvo Fostures

Detection channel

N/A

Call us: 800-304-5357 Email: btinfo@biotium.com

Antibody # prenx	Conjugation	EX/EIII (IIIII)	Laser line	Detection chainer	Dyc i catales
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	

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 3, 2025 at 12:15:02 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/ttf-1-nkx2-1-recombinant-monoclonal-rabbit-antibody-nx2-1-1855r/">https://biotium.com/product/ttf-1-nkx2-1-recombinant-monoclonal-rabbit-antibody-nx2-1-1855r/</a>

N/A

Ev/Em (nm)