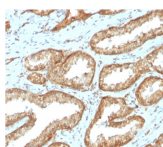


YBX1 / Y-box Binding Protein 1 / YB-1

Monoclonal Mouse Antibody

(YBX1/2430)



Product Description

The specificity of this monoclonal antibody to its intended target was validated by HuProt™ Array, containing more than 19,000 full-length human proteins. Y-box binding protein-1 (YBX1) is the prototypic member of the cold shock protein family that fulfills numerous cellular functions. In the nucleus, YBX1 protein orchestrates transcription of proliferation-related genes, whereas in the cytoplasm it associates with mRNA and directs translation. In human tumor entities, such as breast, lung and prostate cancer, cellular YBX1 expression indicates poor clinical outcome, suggesting that YBX1 is an attractive marker to predict patients' prognosis.

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 2430, Anti-Y-box Binding Protein 1|YB-1|YBX1 (YBX1/2430)

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

Product attributes

Antibody number	#2430
Antibody reactivity (target)	Y-box Binding Protein 1, YB-1, YBX1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	YBX1/2430
Isotype	IgG2b, kappa
Molecular weight	36 kDa
Synonyms	BP-8; CSDA2; CSDB; DBPB; MDR-NF1; NSEP-1; NSEP1; YB-1; YB1
Human gene symbol	YBX1
Entrez gene ID	4904
SwissProt	P67809
Unigene	473583
Immunogen	Recombinant human full-length YBX1 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human
Positive control	MCF-7, HEK293 or PC12 cells. Prostate, Breast or Lung 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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Tumor expression	Breast cancer, Lung cancer, Prostate cancer

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 10, 2025 at 06:15:12 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ybx1-y-box-binding-protein-1-yb-1-monoclonal-mouse-antibody-ybx1-2430/>