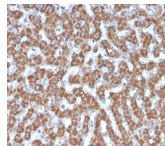


Prohibitin Monoclonal Mouse Antibody (PHB/3226)



Product Description

Recognizes a protein of 30 kDa which is identified as Prohibitin, an evolutionarily conserved protein with homologues found in yeast to man. It is located in the inner membrane of mitochondria. Although prohibitin mRNA and protein expression occurs throughout the cell cycle, maximum levels are detected during the G1/S phase transition and minimum levels are seen in S phase and the G2/mitosis boundary. Prohibitin is located exclusively in the mitochondria with the highest concentration on the inner membrane. Prohibitin is an ideal mitochondrial marker. It shows antiproliferative activity and has been proposed to play a role in normal cell cycle regulation, replicative senescence, cellular immortalization, and tumor suppression. 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.

Product attributes

Antibody number	#3226
Antibody reactivity (target)	Prohibitin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PHB/3226
Isotype	IgG2b, kappa
Molecular weight	~30 kDa
Synonyms	Epididymis luminal protein 215; Epididymis secretory sperm binding protein Li 54e; HEL 215; HEL S 54e; PHB; PHB1; Prohibitin
Human gene symbol	PHB
Entrez gene ID	5245
SwissProt	P35232
Unigene	514303
Immunogen	Recombinant human PHB protein fragment (aa167-261) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Mitochondria
Species reactivity	Human
Positive control	Liver, Kidney, Heart or Endometrium (IHC).
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

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
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, LI-COR Bioscience.