SHBG Monoclonal Mouse Antibody (SHBG/245)



Product Description

Recognizes a protein of 45 kDa, identified as SHBG. It functions as an androgen transport protein, but may also be involved in receptor-mediated processes. Each dimer binds one molecule of steroid. It is specific for 5-alpha-dihydrotestosterone, testosterone, and 17-beta-estradiol. SHBG regulates the plasma metabolic clearance rate of steroid hormones by controlling their plasma concentration. In testis, it is synthesized by the Sertoli cells, secreted into the lumen of the seminiferous tubule and transported to the epididymis.

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 0245, Anti-SHBG (SHBG/245)

Product attributes

Product attributes				
Antibody number	#0245			
Antibody reactivity (target)	SHBG			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	SHBG/245			
Isotype	lgG1			
Molecular weight	45 kDa			
Synonyms	Androgen binding protein (ABP); Sex hormone-binding globulin (SHBG); Sex steroid binding protein (SBP); Sex steroid-binding protein; Testis-specific androgen-binding protein; Testosterone binding beta globulin; Testosterone-Estradiol-Binding Globulin (TEBG)			
Human gene symbol	SHBG			
Entrez gene ID	6462			
SwissProt	P04278			
Unigene	632235			
Immunogen	Recombinant full-length human SHBG protein			
Antibody target cellular localization	Secreted (extracellular)			
Verified antibody applications	IHC (FFPE) (verified)			
Species reactivity	Human			
Antibody application notes	Immunohistology formalin-fixed 1-2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM ris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dillution for a specific application should be determined by user			
Positive control	Liver, Placenta or Testis			
Shipping condition	Room temperature			
Storage Conditions	ore at 2 to 8 °C, Protect fluorescent conjugates from light, ote: 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/m in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL PBS without azide			
Antibody research areas	Endocrinology			
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.			
Cell/tissue expression	Testis			

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

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 4, 2025 at 10:53:10 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-shbg-sex-hormone-binding-globulin-shbg245/