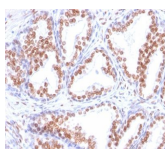


Androgen Receptor Monoclonal Mouse Antibody (DHTR/882)



Product Description

This antibody recognizes a protein of 110 kDa, which is identified as androgen receptor (AR). It reacts with full length, and the newly described A form of the receptor. It does not cross react with estrogen, progesterone, or glucocorticoid receptors. AR functions as a steroid-hormone activated transcription factor. Upon binding the hormone ligand, the receptor dissociates from accessory proteins, translocates into the nucleus, dimerizes, and then stimulates transcription of androgen responsive genes. The expression of AR is reportedly inversely correlated with histologic grade; ie, well differentiated prostate tumors show higher expression than the poorly differentiated tumors. In prostate cancer, AR has been proposed as a marker of hormone-responsiveness and thus it may be useful in identifying patients likely to benefit from anti-androgen therapy. Anti-androgen receptor has been useful clinically in differentiating morpheiform basal cell carcinoma (mBCC) from desmoplastic trichoepithelioma (DTE) in the skin. This MAb is superb for staining of formalin/paraffin tissues.

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 0882, Anti-Androgen Receptor (DHTR/882)

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

Product attributes

Antibody number	#0882
research-areas	Cancer, Endocrinology, Signal transduction
Antibody reactivity (target)	Androgen Receptor
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DHTR/882
Isotype	IgG1, kappa
Molecular weight of antigen	~110 kDa
Synonyms	Androgen Receptor; AR; Dihydrotestosterone Receptor; Nuclear Receptor Subfamily 3 Group C Member 4
Human gene symbol	AR
Entrez gene ID	367
SwissProt	P10275
Unigene	496240
Immunogen	Recombinant full-length human DHTR protein
Antibody target cellular localization	Cytoplasmic, Nucleus
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL. Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Does not react with mouse; others not known. Optimal dilution for a specific application should be determined by user
Positive control	LNcap cells or Prostate 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
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, Endocrinology, Signal transduction
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	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 June 14, 2026 at 01:41:23 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-androgen-receptor-dhtr882/>