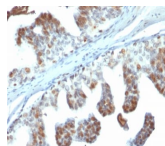


PTEN Monoclonal Mouse Antibody (PTEN/2110)



Product Description

PTEN (phosphatidylinositol-3,4,5-trisphosphate 3-phosphatase) contains a tensin like domain as well as a catalytic domain similar to that of the dual specificity protein tyrosine phosphatases. It was identified as a tumor suppressor that is mutated in a large number of cancers, including sporadic brain, breast, kidney, and prostate cancers. 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 2110, Anti-PTEN (PTEN/2110)**

Product attributes

Antibody number	#2110
Antibody reactivity (target)	PTEN
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PTEN/2110
Isotype	IgG2b, kappa
Molecular weight	55 kDa
Synonyms	BZS; DEC; GLM2; MHAM; Mutated in Multiple Advanced Cancers 1 (MMAC1); MMAC1 phosphatase and tensin homolog deleted on chromosome 10; Phosphatase and Tensin Homolog; Phosphatase and tensin like protein; Phosphatidylinositol 3,4,5-trisphosphate 3-phosphatase and dual-specificity protein phosphatase; PTEN; PTEN1; TEP1
Human gene symbol	PTEN
Entrez gene ID	5278
SwissProt	P60484
Unigene	500466; 729457
Immunogen	Recombinant full-length human PTEN protein
Antibody target cellular localization	Nucleus & cytoplasm
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. Immunohistochemistry (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user
Positive control	MCF-7 or HeLa Cells. Prostate, Breast, Liver, Renal 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
Validated in protein arrays	Monospecific
Antibody research areas	Signal transduction, Tumor suppressors

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

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.