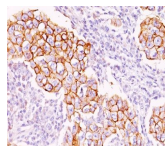


Phosphotyrosine Monoclonal Mouse Antibody (PY20)



Product Description

Protein phosphorylation is a fundamental event in the regulation of a large number of intracellular processes. Phosphorylation of specific tyrosine residues is the result of activation or stimulation of their respective protein tyrosine kinases. The phosphorylated proteins can be auto-phosphorylated kinases or certain cellular protein substrates. Tyrosine-phosphorylated proteins are involved in signal transduction and in the regulation of cell proliferation. Antibody to phosphotyrosine provides an excellent tool for the detection, characterization, and purification of phosphotyrosine containing proteins. This MAb shows no cross-reaction with other phosphoamino acids and is superb for multiple applications including 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 0232, Anti-Phosphotyrosine (PY20)**

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Biol Chem (1983) 264(34): 20163-20166. (IP)
2. J Immunological Meth (1988) 109: 277-285. (ELISA)
3. Mol Endocrinol (1998) 12(4): 536-543. (WB)
4. Fert Stert (1999) 71(5): P919-923. (IF)

Product attributes

Antibody number	#0232
Antibody reactivity (target)	Phosphotyrosine
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PY20
Isotype	IgG2b
Molecular weight	Depends upon the phosphorylated target
Synonyms	P-Tyr
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
UniGene	Not Applicable
Immunogen	Phosphotyrosine conjugated to KLH
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	All species
Expected antibody applications	ELISA (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL Immunohistology formalin-fixed 1-2 ug/mL. No pretreatment is required for staining of formalin-fixed, paraffin-embedded tissues. Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1 ug/mL, Immunoprecipitation 1-2 ug/500 ug protein lysate, Optimal dilution for a specific application should be determined by user
Positive control	MCF-7, MDA-231, T47-D cells or breast 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	Signal transduction

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.

This datasheet was generated on May 18, 2024 at 05:58:51 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-phosphotyrosine-py20/>