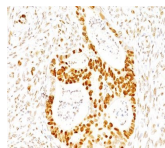


p27 / KIP1 Monoclonal Mouse Antibody (SX53G8)



Product Description

This MAb recognizes a 27 kDa protein, identified as the p27Kip1, a cell cycle regulatory mitotic inhibitor. It is highly specific and shows no cross-reaction with other related mitotic inhibitors. In Western blotting of cell lysates from 7 human breast cancer cell lines (ZR75-1, ZR75-30, MCF-7, MDAMB453, T47D, CAL51, 734B), the antibody labels a single band corresponding to p27Kip1. It functions as a negative regulator of G1 progression and has been proposed to function as a possible mediator of TGF- β -induced G1 arrest. p27Kip1 is a candidate tumor suppressor gene. Reportedly, low p27 expression has been associated with unfavorable prognosis in renal cell carcinoma, colon carcinoma, breast carcinomas, non-small-cell lung carcinoma, hepatocellular carcinoma, multiple myeloma, and lymph node metastases in papillary carcinoma of the thyroid, as well as a more aggressive phenotype in carcinoma of the cervix. Primary antibodies are available purified, or with a selection of fluorescent CF Φ dyes and other labels. CF Φ dyes offer exceptional brightness and photostability. See the [CF \$\Phi\$ Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF Φ 405S and CF Φ 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 Φ 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 0669, Anti-KIP1 (SX53G8)**

References

Note: References for this clone sold by other suppliers may be listed for expected applications. 1. PNAS USA (1997) 94:6380-6385. (western; immunoprecipitation; IHC, FFPE) 2. Blood (2005) 105(9): 3691. (western) 3. Anticancer Res (2017) 37: 2407-2415. (IHC, FFPE; IF, western)

Product attributes

Antibody number	#0669
Antibody reactivity (target)	KIP1, p27
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SX53G8
Isotype	IgG1, kappa
Molecular weight	25-26 kDa
Synonyms	CDKN1B, CDKN4, Cyclin Dependent Kinase Inhibitor 1B, Cyclin-dependent kinase inhibitor p27 Kip1, KIP1, MEN1B, MEN4
Human gene symbol	CDKN1B
Entrez gene ID	1027
SwissProt	P46527
Unigene	238990
Immunogen	Purified GST-p27 fusion protein of human origin
Antibody target cellular localization	Nucleus
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified)
Species reactivity	Human, Monkey, Mouse, Rat
Expected antibody applications	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: 0.5-1 μ g/mL. Immunohistochemistry formalin-fixed 0.25-0.5 μ g/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 μ g/million cells/0.1 mL. Western blotting 0.5-1 μ g/mL. Optimal dilution for a specific application should be determined by user
Positive control	ZR75, T47D, SK-BR-3, MDA-MB-231, HeLa or MCF7 cells. Tonsil, Breast, Cervical or Colon Carcinoma.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 $^{\circ}$ C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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	Cell cycle, Tumor suppressors

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