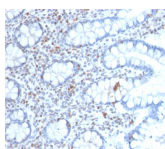


BMI1 Monoclonal Mouse Antibody (BMI1/2689)



Product Description

The B cell-specific moloney murine leukemia virus integration site 1 (Bmi-1) is a transcriptional receptor of the polycomb gene family involved in several cellular processes including cell-cycle regulation, apoptosis, and maintenance of adult and neoplastic stem cells by providing self-renewal capacity. Further, Bmi-1 expression has been associated with malignant transformation, tumor progression, metastatic disease, and poor prognosis in human malignancies.

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 2689, Anti-BMI1 (BMI1/2689)

Product attributes

Antibody number	#2689
Antibody reactivity (target)	BMI1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BMI1/2689
Isotype	IgG1, kappa
Molecular weight	~41 kDa
Synonyms	BMI1; Polycomb complex protein BMI1; B lymphoma Mo MLV insertion region 1 homolog; BMI1 polycomb ring finger oncogene; Polycomb group ring finger 4; RING finger protein 51
Human gene symbol	BMI1
Entrez gene ID	648
SwissProt	P35226
Unigene	380403
Immunogen	Recombinant fragment of human BMI1 protein (around aa 142-326) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoplasmic, Nucleus
Species reactivity	Human
Positive control	293T, NIH3T3, HeLa or Jurkat cells. Colon 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
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 array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody research areas	Cancer, Developmental biology
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.

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 September 10, 2025 at 10:36:33 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bmi1-monoclonal-mouse-antibody-bmi1-2689/>