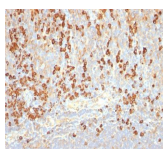


# Plasma Cell Marker Monoclonal Mouse Antibody (LIV3G11, same as 7B18)



## Product Description

This antibody recognizes an intra-cytoplasmic antigen, which shows a very high degree of specificity for plasma cells. This antigen is present in normal as well as neoplastic plasma cells. Plasma cells, which are large lymphocytes derived from an antigen-specific B cell, secrete antibodies and are responsible for humoral immunity. Plasma cells differentiate from B cells upon stimulation by CD4 lymphocytes. The B cell acts as an antigen-presenting cell (APC), consuming an offending pathogen, which is taken up by the B cell by phagocytosis and broken down within proteosomes. Plasma cells contain basophilic cytoplasm; their nucleus contains heterochromatin organized in a characteristic cartwheel arrangement. This MAb superbly recognizes normal and neoplastic plasma cells in routine formalin-fixed, paraffin-embedded tissue sections. It is of potential value in identifying myeloma or plasmacytoma in bone marrow or other tissues. It also helps differentiate lympho-plasmacytoid lymphoma from lymphocytic and follicular lymphoma. Note that this MAb is not suitable for staining frozen 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0047, Anti-Plasma Cell Marker (LIV3G11; same as 7B18)**

## Product attributes

Antibody number	#0047
Antibody reactivity (target)	Plasma Cell Marker
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LIV3G11; same as 7B18
Isotype	IgG2a
Molecular weight	Not Known
Synonyms	Plasma B cell, Plasmocytes
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Pancreatic cancer-related mucin
Antibody target cellular localization	Cytoplasmic
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Expected antibody applications	IF (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. Does not react with rat; others not known. Immunohistology formalin-fixed 0.1-0.2 ug/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. Optimal dilution for a specific application should be determined by user
Positive control	Tonsil or lymph node
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	Immunology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	B-cells, Lymphocytes
Tumor expression	Leukemia/lymphoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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.

## References

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

Front Immunol (2017) 8:563. (IF, FFPE)

This datasheet was generated on January 10, 2025 at 01:26:20 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-plasma-cell-marker-liv3g11-7b18/>