MCM7 Recombinant Monoclonal Rabbit Antibody (MCM7/2756R)



Product Description

MCM7 is one of the highly conserved mini-chromosome maintenance proteins (MCM) that is essential for the initiation of eukaryotic genome replication. The hexameric protein complex formed by the MCM proteins is a key component of the pre-replication complex and may be involved in the formation of replication forks and in the recruitment of other DNA replication related proteins. The MCM complex consisting of this protein and MCM2, 4 and 6 proteins possesses DNA helicase activity, and may act as a DNA unwinding enzyme. Cyclin D1-dependent kinase, CDK4, is found to associate with this protein, and may regulate the binding of this protein with the tumor suppressor protein RB1/RB.

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 2756, Anti-MCM7 (MCM7/2756R)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#2756			
Antibody reactivity (target)	MCM7			
Antibody type	Primary			
Host species	Rabbit			
Clonality	Recombinant Monoclonal			
Clone	MCM7/2756R			
Isotype	IgG			
Molecular weight	88 kDa			
Synonyms	CDC47; DNA replication licensing factor MCM7; MCM7 mini chromosome maintenance deficient 7; Mini chromosome Maintenance 7; Mini chromosome maintenance protein 7; P1.1-MCM3; P1CDC47; P85MCM; PNAS146			
Human gene symbol	MCM7			
Entrez gene ID	4176			
SwissProt	P33993			
Unigene	438720			
Immunogen	Recombinant fragment of human MCM7 (aa195-319)			
•				
Verified antibody applications	IHC (FFPE) (verified), WB (verified)			
	· , ,			
applications Antibody target cellular	IHC (FFPE) (verified), WB (verified)			
applications Antibody target cellular localization	IHC (FFPE) (verified), WB (verified) Nucleus			
applications Antibody target cellular localization Species reactivity	IHC (FFPE) (verified), WB (verified) Nucleus Human			
applications Antibody target cellular localization Species reactivity Positive control	IHC (FFPE) (verified), WB (verified) Nucleus Human Raji or Ramos cells. Tonsil			
applications Antibody target cellular localization Species reactivity Positive control Shipping condition	IHC (FFPE) (verified), WB (verified) Nucleus Human Raji or Ramos cells. Tonsil Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions	IHC (FFPE) (verified), WB (verified) Nucleus Human Raji or Ramos cells. Tonsil Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	IHC (FFPE) (verified), WB (verified) Nucleus Human Raji or Ramos cells. Tonsil Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 m			

Chinese hamster ovary cells. Inquire for the specific lot.

Email: techsupport@biotium.com

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 October 28, 2025 at 07:35:21 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mcm7-recombinant-monoclonal-rabbit-antibody-mcm7-2756r/