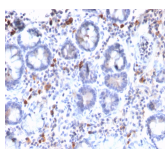


Mucin 3 / MUC 3 Recombinant Monoclonal Rabbit Antibody (MUC3/2992R)



Product Description

Recognizes a protein of HMW, identified as mucin 3 glycoprotein (MUC3). Its epitope localizes between aa SITTTTE. This MAb shows no cross-reaction with human milk fat globule membranes, MUC1, or MUC2. MUC3 is distributed in colon and rectum, and is also present to a lesser extent in breast, lung and salivary gland tissues. The Mucins are a family of highly glycosylated, secreted proteins with a basic structure consisting of a variable number of tandem repeats (VNTRs) encoded by 60 base pairs (Mucin 1), 69 base pairs (Mucin 2) and 51 base pairs (Mucin 3). The number of repeats is highly polymorphic and varies among different alleles. Mucin 1 proteins are expressed as type I membrane proteins in addition to secreted forms. Mucin 1 is aberrantly expressed in epithelial tumors including breast carcinomas. Mucin 2 coats the epithelia of the intestines and airways and is associated with colonic tumors. Mucin 3 is a major component of various mucus gels and is broadly expressed in normal and tumor cells.

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 2992, Anti-MUC3 (MUC3/2992R)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2992
Antibody reactivity (target)	MUC3
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	MUC3/2992R
Isotype	IgG
Molecular weight	1100 kDa
Synonyms	Intestinal mucin-3A; intestinal mucin-like; MUC3; MUC3A; MUC3B; mucin 3; intestinal; mucin 3A; cell surface associated; mucin 3A; intestinal; mucin 3B; cell surface associated
Human gene symbol	MUC3A
Entrez gene ID	4584 & 57876
SwissProt	Q02505 & Q9H195
Unigene	744422 & 744530
Immunogen	A synthetic peptide of 35 amino acids, SIB35 (C-HSTPSFTSSITTTETTSHTSPSTSSITTTETTS), which contains two of the MUC3 tandem repeats, coupled to KLH.
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	LS174T cells. Human small intestine (IHC).
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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.
Cell/tissue expression	Epithelial cells, Intestine
Antibody research areas	Cancer, Mucins

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 December 11, 2025 at 03:19:00 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mucin-3-muc-3-recombinant-monoclonal-rabbit-antibody-muc3-2992r/>