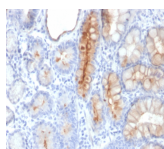


# Mucin 4 / MUC4 Monoclonal Mouse Antibody (MUC4/3084)



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

## Product Description

The major constituents of mucus, the viscous secretion that covers epithelial surfaces such as those in the trachea, colon, and cervix, are highly glycosylated proteins called mucins. These glycoproteins play important roles in the protection of the epithelial cells and have been implicated in epithelial renewal and differentiation. This gene encodes an integral membrane glycoprotein found on the cell surface, although secreted isoforms may exist. MUC-4 transcripts have been detected in normal respiratory epithelium and lung. MUC-4 is a very specific (100%) and sensitive (90%) marker of lung adenocarcinomas and is negative for mesotheliomas. Reportedly, MUC-4 expression in invasive ductal carcinoma of the pancreas is an independent factor for poor prognosis.

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 3084, Anti-MUC4 (MUC4/3084)**

## Product attributes

Antibody number	#3084
Antibody reactivity (target)	MUC4
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MUC4/3084
Isotype	IgG1, kappa
Molecular weight of antigen	235 kDa
Synonyms	Ascites sialoglycoprotein 1 / 2; ASGP; ASGP-1; ASGP-2; MUC-4; MUC4; Mucin 4 cell surface associated; Mucin 4 tracheobronchial; Mucin-4 beta chain; Pancreatic adenocarcinoma mucin; Testis mucin; Tracheobronchial mucin
Human gene symbol	MUC4
Entrez gene ID	4585
SwissProt	Q99102
Unigene	369646
Immunogen	Recombinant fragment (around aa 1730-1864) of human MUC4 protein (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Positive control	Stomach, Colon or esophagus.
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
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	Epithelial cells
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	<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.

This datasheet was generated on June 19, 2026 at 08:53:24 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mucin-4-muc4-monoclonal-mouse-antibody-muc4-3084/>