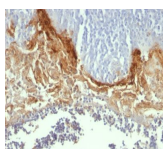


Involucrin Monoclonal Mouse Antibody (SY5)



Product Description

This antibody recognizes a protein of 66 kDa-170 kDa, identified as involucrin. In Western blotting of cultured human keratinocytes, this MAb reacts with a 120 kDa protein. It stains the involucrin in a variety of sizes: 170 kDa in MCF-7 cells, a doublet of ~115 kDa and 150 kDa in gorilla and owl monkey, 66 kDa in dog, and a doublet of 105 kDa in pig. Its epitope maps between codon 421-568 of human involucrin. Involucrin is expressed in a range of stratified squamous epithelia, including the cornea, which lacks a distinct cornified layer. In normal epidermis, it is first expressed in the upper spinous layers, and in keratinocyte cultures, all cells that have left the basal layer express it. Involucrin expression is altered in pathological conditions: in psoriasis and other benign epidermal hyperplasias, involucrin expression begins closer to the basal layer than normal; expression is abnormal in squamous cell carcinomas and premalignant lesions, and is reduced in severe dysplasias of the larynx and cervix.

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 0678, Anti-Involucrin (SY5)

Product attributes

Antibody number	#0678
Antibody reactivity (target)	Involucrin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SY5
Isotype	IgG1, kappa
Molecular weight of antigen	66-170 kDa
Synonyms	IVL
Human gene symbol	IVL
Entrez gene ID	3713
SwissProt	P07476
Unigene	516439
Immunogen	Purified involucrin from human keratinocytes
Antibody target cellular localization	Cytoplasmic
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Dog, Gorilla, Human, Monkey, Owl, Pig
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology formalin-fixed 0.1-0.2 ug/mL. Staining of formalin-fixed tissues is enhanced by digestion with trypsin or Protease XXV at 1mg/mL PBS for 5 min at 37C, enzyme digestion is better than citrate buffer based epitope unmasking. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Immunofluorescence. Does not react with mouse; others not known. Optimal dilution for a specific application should be determined by user
Positive control	MCF-7 cells. Localized to upper spinous and granular layers in normal skin.
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	Cancer
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.
Tumor expression	Squamous cell carcinoma

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 June 15, 2026 at 06:40:51 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-involucrin-sy5/>