

# MUC1 Recombinant monoclonal antibody (SM3) (Mouse IgG1 $\lambda$ )

This antibody is covered by our [Worry-Free Guarantee](#).

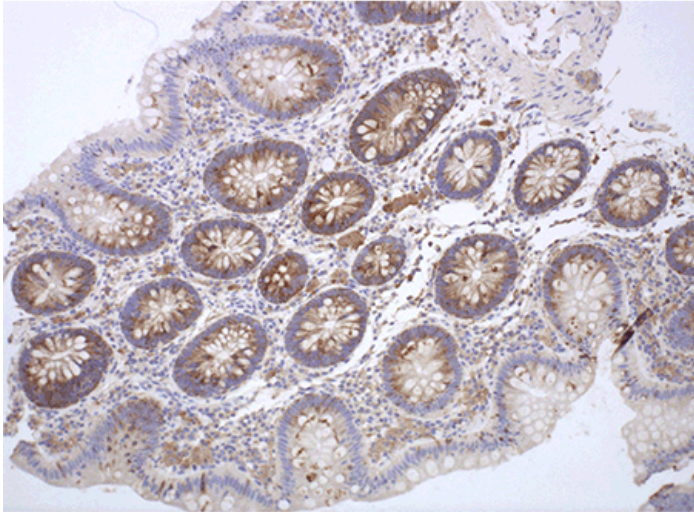
## Ordering Information

[Order Online »](#)

ENZ-ABS645-0200	200 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded colon tissue stained using Enzo's MUC1 antibody (Prod. No. ENZ-ABS645). IHC staining was performed on an automated Leica BOND III.

# Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Mucin
Application	ELISA, Flow Cytometry, IHC (FS), IHC (PS)
Clone	SM3
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Hydrogen fluoride deglycosylated milk mucin.
Isotype	IgG1λ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Recognizes the under-glycosylated form of human MUC1, a marker of tumors.
UniProt ID	P15941
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)