

MUC1 Recombinant monoclonal antibody (SM3) (Mouse IgG1 λ)

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

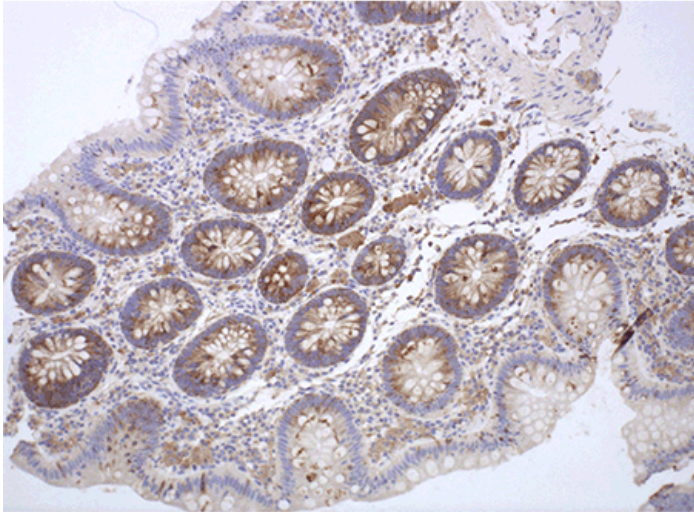
Ordering Information

[Order Online »](#)

ENZ-ABS645-0200	200 μ 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	IgG1 λ
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 [Worry-Free Guarantee](#).



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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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