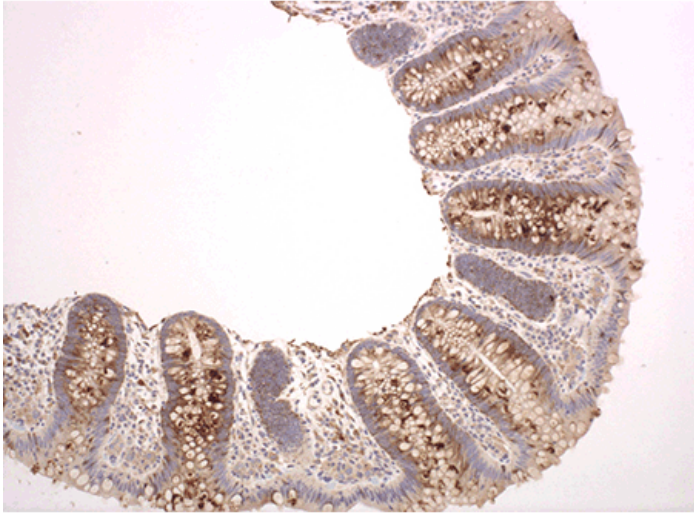


MUC1 Recombinant monoclonal antibody (HMFG2) (Rabbit IgG λ)

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

Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded colon tissue stained using Enzo's MUC1 antibody (Prod. No. ENZ-ABS644). 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, IHC
Clone	HMFG2
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	High-MW glycoprotein of the human milk fat globule.
Isotype	IgG λ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original Mouse IgG1 format, for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P15941
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



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)