

MUC1 Recombinant monoclonal antibody (rMUC1/960)

Mucin-1/MUC1 protein is encoded by the MUC1 gene in humans. This gene encodes a membrane-bound protein that is a member of the mucin family. Family members are O-glycosylated proteins that play an essential role in forming protective mucous barriers on epithelial surfaces. MUC1 is a transmembrane protein normally expressed on the apical borders of secretory epithelial cells. Overexpression is often associated with colon, breast, ovarian, lung and pancreatic cancers.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Mucin-1, EMA, Tumor-associated epithelial membrane antigen, CD227
Application	Flow Cytometry, IF, IHC (PS)
Clone	rMUC1/960
Formulation	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Gene/Protein Identifier	4582 (Gene ID)
Host	Mouse
Immunogen	Human milk fat globule membranes.
Isotype	IgG2bk
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1 µg/10 ⁶ cells in 0.1 mL)Immunofluorescence (1-2 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (1-2 µg/mL for 30 minutes at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
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)