MUC1 Recombinant monoclonal antibody (HMFG2) (Mouse IgG1λ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS643-0200

200µg

Manuals, SDS & CofA

View Online »

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 Mouse

Immunogen High-MW glycoprotein of the human milk fat globule.

lsotype IgG1λ

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P15941

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com