

MUC1 Recombinant monoclonal antibody (HMFG2) (Mouse IgG1 λ)

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

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.
Isotype	IgG1 λ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P15941
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee