Tumor associated glycoprotein TAG-72 Recombinant monoclonal antibody

(umurasociate) glycopydar 20 9 dlycopythin on the surface of many cancer cells, including ovary, breast, colon, and pancreatic cells.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS567-0200

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

Application Flow Cytometry, IHC

Clone B72.3 (Satumomab)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Human tumor associated glycoprotein 72.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and

techniques.

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

