HLA-G (human) monoclonal antibody (MEM-G/1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

Ordering Information

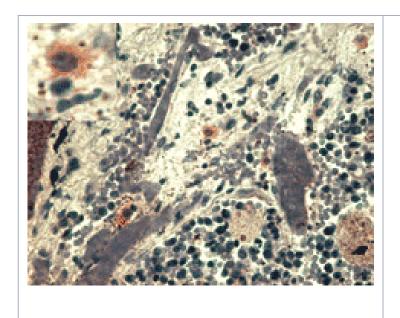
Order Online »

ALX-805-704-C100

100µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Handling Keep at +4°C. Do not freeze.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Clone MEM-G/1

Formulation Liquid. In PBS, pH 7.4, containing 15mM sodium azide.

Host Mouse

Immunogen Human recombinant HLA-G denatured heavy chain.

lgG1

Purity Detail Protein A affinity chromatography.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, stain by standard

ABC technique, pretraet by heat retrieval in 0.01 citrate buffer (4 x 2 minutes in microwave oven); recommended starting dilution is 1:60-1:100; incubate 1 hour at room

temperature; positive tissue: Human placenta –

extravillous cytotrophoblast)Suggested dilutions/conditions may not be available for all applications.Optimal conditions

must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes denatured HLA-G heavy chain.

UniProt ID P17693

Last modified: May 29, 2024



info-

eu@enzolifesciences.com