IL-12 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

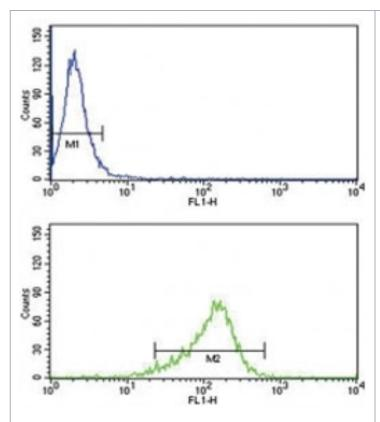
Order Online »

ENZ-ABS895-0400

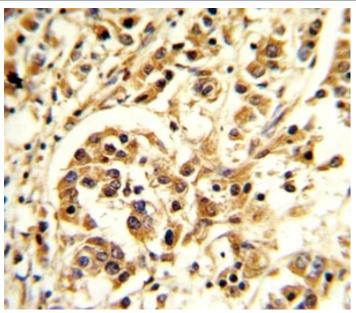
0.4ml

Manuals, SDS & CofA

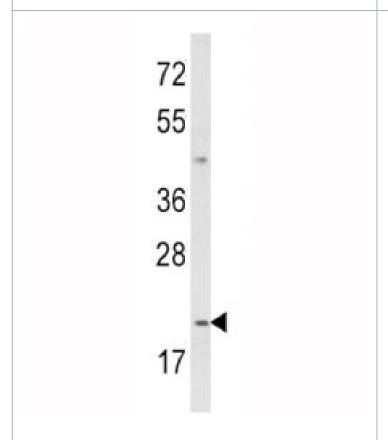
View Online »



IL-12 antibody flow cytometry analysis of MDA-MB231 cells (green) compared to a negative control (blue). FITC-conjugated goat-anti-rabbit secondary Ab was used for the analysis.



IHC analysis of FFPE human breast carcinoma stained with IL-12 antibody



Western blot analysis of IL-12 antibody and MDA-MB231 lysate. Predicted molecular weight ~ 24 kDa.

Handling & Storage

Handling Aliquot the IL-12 antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-12

Application ELISA, Flow Cytometry, IHC, WB

Formulation Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.

Host Rabbit

Immunogen Human IL-12 (aa 168-195).

Isotype Ig

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin

sections) (1:10-1:50)Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P29459

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

