

IFNA1 antibody (mAb)

Catalog Nos: 61603, 61604

RRID: AB_2793697

Isotype: IgG2a

Application(s): ELISA

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 0.82 µg/µl

Molecular Weight: 19 kDa

Background: IFNA1 (Interferon, Alpha 1) is produced by macrophages, IFN-alpha have antiviral activities. Interferon stimulates the production of two enzymes: a protein kinase and an oligoadenylate synthetase.

Immunogen: Human IFNA1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

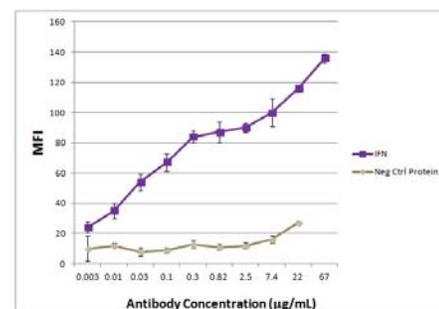
Application Notes:

Applications Validated by Active Motif:

ELISA: Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



IFNA1 antibody (mAb) tested by ELISA.

Recombinant IFNA1 protein or a negative control protein was conjugated to Luminex® microspheres and then incubated in the same well with a serial dilution of IFNA1 antibody (mAb). Biotinylated anti-mouse IgG followed by Streptavidin-PE was used to detect the bound IFNA1 antibody. Mean Fluorescence Intensity (MFI) for each protein was measured in a Luminex® 200™.