

GATA-4 antibody (pAb)

Catalog Nos: 39893, 39894

RRID: AB_2615046

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.66 μ g/ μ l

Molecular Weight: 55 kDa

Background: GATA-4 is a zinc finger protein and part of the GATA family of sequence specific DNA binding transcription factors that recognize a consensus DNA sequence containing GATA. GATA-4 is involved in the development of several mesodermal and endodermal structures. GATA-4 mutant mice exhibit severe heart malformations. Defects in GATA-4 are the cause of atrial septal defect type 2. ASD2 is a congenital heart malformation characterized by incomplete closure of the wall between the atria resulting in blood flow from the left to the right atria. GATA-4 also plays a role in regulating gene expression in cardiac and smooth muscle cell types.

Immunogen: This GATA-4 antibody was raised against a peptide derived from the C-terminus of human GATA-4.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

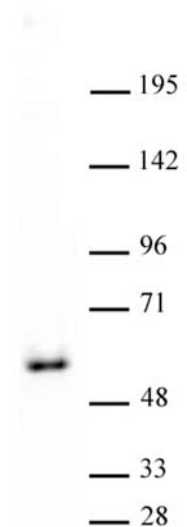
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 dilution

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.



GATA-4 antibody (pAb) tested by Western blot.

Western Blot: Nuclear extract (20 μ g) of F9 cells (retinoic acid treated mouse teratocarcinoma F9 stem cell line, Active Motif Catalog no. 36113) probed with GATA-4 antibody at a dilution of 1:2,000.