

GATA-6 antibody (pAb)

Catalog Nos: 61063, 61064

RRID: AB_2793493

Isotype: IgG

Application(s): ICC, IF, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 62 kDa

Background: GATA-6 (GATA binding factor 6) is a Zn-finger DNA binding transcription factor that is required for the expression of genes regulating formation of the primitive endoderm lineage. GATA-6, along with GATA-4, is a marker of extraembryonic endoderm.

Immunogen: This GATA-6 antibody was raised against a peptide within the N-terminus of human GATA-6.

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:

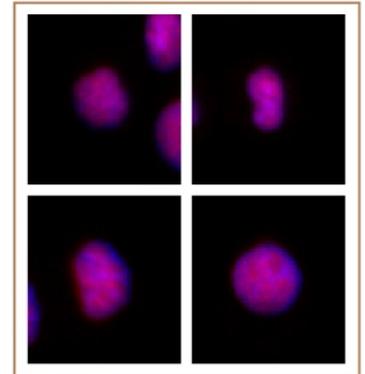
Applications Validated by Active Motif:

ICC/IF: 1:2,000 dilution

WB: 1:500 dilution

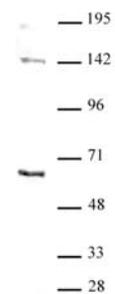
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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-6 antibody (pAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with GATA-6 antibody at a 1:2,000 dilution.



GATA-6 antibody (pAb) tested by Western blot.

Nuclear extract (20 μ g) of SW48 cells probed with GATA-6 antibody at a dilution of 1:500.