

Anti-GATA1

Rabbit Polyclonal Antibody

Catalog # G19-63R

Lot # R157-5

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the GATA1 protein

Cross Reactivity

- Western blot human and mouse GATA1

GATA1 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 308-312 of human GATA1

Formulation

PBS, pH 7.4, 0.02% NaN₃, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

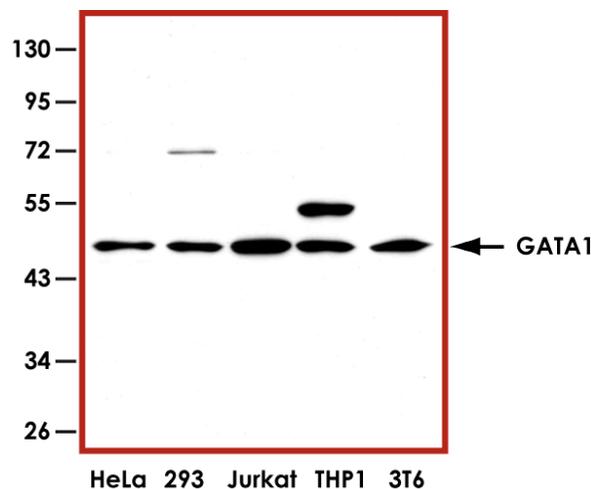
GATA1 is a member of the GATA family of transcription factors and plays an important role in erythroid development by regulating the switch of fetal hemoglobin to adult hemoglobin. FOG, a zinc finger protein, interacts with the amino (N) finger of GATA1 and this interaction is essential for the function of GATA1 in erythroid differentiation (1). Caspase-mediated cleavage of GATA1 is an important negative control mechanism in erythropoiesis. Mutations in GATA1 have been associated with X-linked dyserythropoietic anemia and thrombocytopenia.

References

1. Crispino, J D. et al: Use of altered specificity mutants to probe a specific protein-protein interaction in differentiation: the GATA-1:FOG complex. *Molec. Cell* 3: 219-228, 1999.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Sample Data



Representative Western blot with Anti-GATA1 (1:1000) using 20 µg of HeLa, 293, Jurkat, THP1 and 3T6 cell lysate.

Anti-GATA1

Rabbit Polyclonal Antibody

Catalog Number

G19-63R

Specific Lot Number

R157-5

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.