

## GATA-1 antibody (pAb)

---

**Catalog Nos:** 61535, 61536

**RRID:** AB\_2793673

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Concentration:** 1 µg/µl

**Molecular Weight:** 47 kDa

---

**Background:** **GATA-1** (Globin transcription factor) is a Transcriptional activator which probably serves as a general switch factor for erythroid development. It binds to DNA sites with the consensus sequence [AT]GATA[AG] within regulatory regions of globin genes and of other genes expressed in erythroid cells.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human GATA-1.

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

**Application Notes:**

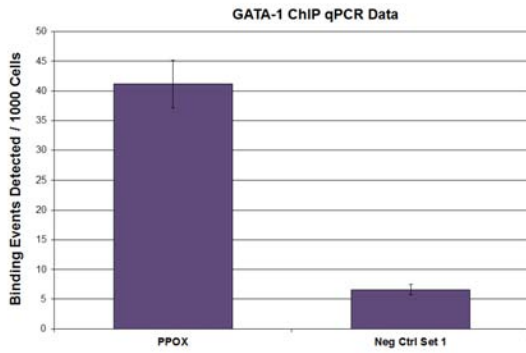
Applications Validated by Active Motif:

ChIP-Seq: 4 µg per ChIP

ChIP: 10 µg per ChIP

**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-1 (pAb) tested by ChIP.

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 15 µg of K562 cells and 10 µg of GATA-1 antibody. ChIP DNA was used in qPCR with the gene-specific primer for PPOX (a gene involved in heme biosynthesis that is regulated by GATA-1) or negative control primer as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.

#### GATA-1 antibody (pAb) tested by ChIP-Seq

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 µg of K562 cells and 4 µg of GATA-1 antibody. ChIP DNA was sequenced on the Illumina NextSeq and 12 million sequence tags were mapped to identify GATA-1 binding sites on chromosome 1.

