

YY1 antibody (pAb)

Catalog No: 39071

Isotype: IgG

Application(s): ChIP, WB

Reactivity: Human, Monkey, Mouse, Rat

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.3 µg/µl

Volume: 200 µl

Molecular Weight: 63 kDa

Background: YY1 (Yin Yang 1) is a transcription factor that belongs to the GLI-Kruppel class of zinc finger transcription factors and is ubiquitously expressed. YY1 is known to have a fundamental role in normal biologic processes such as embryogenesis, differentiation, replication and cellular proliferation. YY1 exerts its effects on genes involved in these processes via its ability to initiate, activate or repress transcription depending upon the context in which it binds. Since expression and function of YY1 are known to be intimately associated with progression through phases of the cell cycle, the physiologic significance of YY1 activity has been applied to models of tumor biology. It can repress and activate a number of gene promoters and direct histone deacetylases (HDACs) and histone acetyltransferases (HATs).

Immunogen: This YY1 antibody was raised against a synthetic peptide corresponding to amino acid residues 109-123 of human YY1.

Buffer: PBS containing 0.2% gelatin, 0.01% thimersol and 0.1% sodium azide. Thimersol and sodium azide are highly toxic.

Application Notes:

Validated Applications:

ChIP: 10-15 µl per ChIP

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

Published Applications:

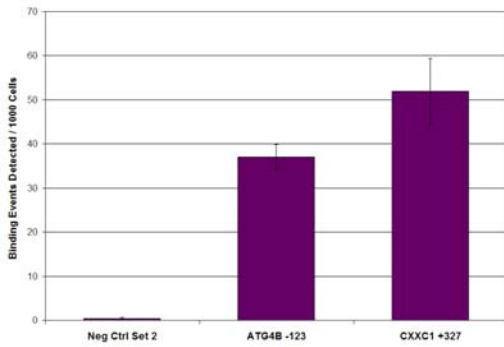
WB

See references for more information. Individual optimization may be required.

For optimal results, primary antibody incubations should be performed at 4°C. Individual optimization may be required.

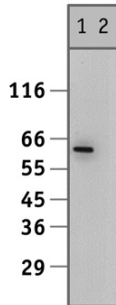
Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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



YY1 antibody tested by ChIP.

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from human lung melanoma cell line MALME-3M cells and 13 µl of YY1 antibody. ChIP DNA was used in qPCR with the control primer pairs or gene-specific primer pairs 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.



YY1 antibody tested by Western blot.

Detection of YY1 by Western blot. The analysis was performed using Raji nuclear extract and YY1 antibody at a 1:1,000 dilution (lane 1). The signal corresponding to YY1 is eliminated by incubating the reaction mixture together with the immunizing peptide (lane 2).