

AML-3/Runx2 antibody (pAb)

Catalog No: 39302

RRID: AB_2793223

Isotype: Serum

Application(s): EMSA, ICC

Reactivity: Human

Volume: 200 µl

Purification: None

Host: Rabbit

Molecular Weight: 56 kDa

Background: AML-3 (acute myeloid leukemia 3 protein) also known as Runx2 (runt-related transcription factor 2) is a multifunctional transcription factor that can either activate or repress transcription. The alpha subunit can bind DNA as a monomer, or DNA binding can be increased by heterodimerization. AML-3 is necessary for osteoblastic differentiation and skeletal morphogenesis.

Immunogen: This AML-3/Runx2 antibody was raised against a synthetic peptide corresponding to amino acid residues 333-348 of human AML-3/Runx2.

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

Application Notes:

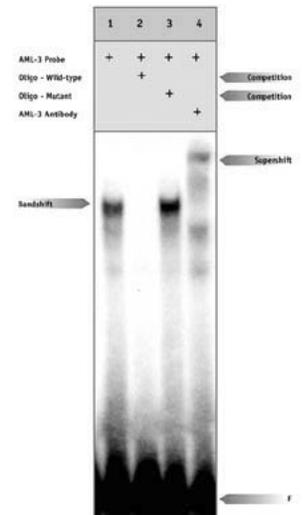
Applications Validated by Active Motif:

EMSA: 2 - 6 µl per assay

The addition of Tris-glycine electrophoresis buffer for supershift assays is recommended.

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.



AML-3/Runx2 pAb tested by supershift.

Detection of AML-3/Runx2 by EMSA analysis. AML-3/Runx2 pAb was used in a supershift assay (EMSA) with Saos-2 nuclear extract (lane 4).