

AML-2/Runx3 antibody (pAb)

Catalog No: 39301

RRID: AB_2793222

Isotype: Serum

Application(s): ChIP, EMSA, ICC, IHC, WB

Reactivity: Human

Volume: 200 μ l

Purification: None

Host: Rabbit

Molecular Weight: 44 kDa

Background: AML-2 (acute myeloid leukemia 2 protein) also known as Runx3 (Runt-related transcription factor 3) is a member of the DNA binding alpha subunit of the heterodimeric transcription factor complex PEBP2/CBF. AML-2 can either activate or repress transcription, and it also interacts with other transcription factors. AML-2 expression is regulated by AML-1/Runx1.

Immunogen: This AML-2 antibody was raised against a full-length protein corresponding to human AML-2/Runx3.

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

Application Notes:

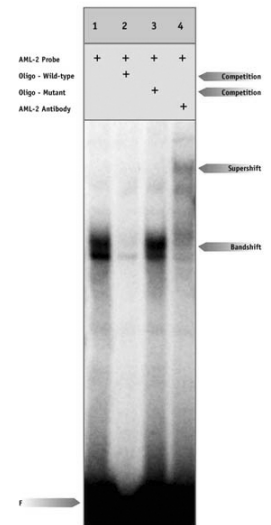
Applications Validated by Active Motif:

EMSA: 4 - 6 μ l 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-2/Runx3 antibody tested by supershift.
Detection of AML-2/Runx3 by EMSA analysis. AML-2/Runx3 antibody was used in a supershift assay (EMSA) with Raji nuclear extract (lane 4).