

Anti-AXL

Rabbit Polyclonal Antibody

Catalog # A34-63R

Lot # Z2014-29

Cited Applications

ELISA, WB

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

Specificity

Recognizes the AXL protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Peptide at the N-terminal residues of human AXL protein.

Formulation

PBS, pH 7.4 with 0.02% Sodium Azide

Stability

1yr at -20°C from date of shipment

Scientific Background

AXL is a member of the receptor tyrosine kinase family which has oncogenic potential and is implicated in human myeloid leukemia (1). AXL is a member of a complex signaling network that is involved in the control of cell proliferation and differentiation. Overexpression of AXL cDNA in NIH 3T3 cells induces neoplastic transformation of these cells with the concomitant appearance of a 140kDa AXL tyrosine-phosphorylated protein (2). Expression of AXL cDNA in the baculovirus system results in the expression of the appropriate recombinant protein that is recognized by anti-phosphotyrosine antibodies, confirming that the AXL protein is tyrosine phosphorylated.

References

1. O'Bryan, J.P. et al: Axl, a transforming gene isolated from primary human myeloid leukemia cells, encodes a novel receptor tyrosine kinase. *Mol Cell Biol.* 1991 Oct;11(10):5016-31.
2. Janssen, J.W. et al.; A novel putative tyrosine kinase receptor with oncogenic potential. *Oncogene.* 1991 Nov;6(11):2113-20.

Anti-AXL

Rabbit Polyclonal Antibody

Catalog # A34-63R

Lot # Z2014-29

Purification Affinity chromatography

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.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.