

Anti-Src

Rabbit Polyclonal Antibody

Catalog # S19-63BR

Lot # Z2014-44

Cited Applications

WB, IHC

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

Specificity

Recognizes the Src protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antiserum was produced against synthesized non-phosphopeptide derived from human Src around the phosphorylation site of tyrosine 529 (P-Q-YP-Q-P).

Formulation

Phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Stability

1yr at -20°C from date of shipment

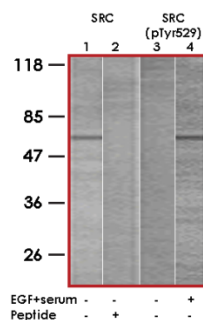
Scientific Background

Src family belongs to non-receptor tyrosine kinases. Src was originally identified as a transforming protein of the Rous sarcoma virus (RSV) that had enzymatic ability to phosphorylate tyrosine in protein substrates (1). Src is overexpressed and activated in a large number of human malignancies and has been linked to the development of cancer and progression to distant metastases (2). In addition to increasing cell proliferation, a key role of Src in cancer seems to be the ability to promote invasion and motility, functions that might contribute to tumour progression.

References

- Collett, M S. et al: Protein kinase activity associated with the avian sarcoma virus src gene product. Proc Natl Acad Sci U S A. 1978 Apr;75(4):2021-4.
- Jacobs, C. et al: Expression of pp60c-src protein kinase in adult and fetal human tissue: high activities in some sarcomas and mammary carcinomas. Cancer Res. 1983 Apr;43(4):1696-702.

Sample Data



Western Blot analysis of extracts from 293 cells using Anti-Src (Lane 1 and 2) and Anti-phospho-Src (Tyr529) (Lane 3 and 4).

Anti-Src

Rabbit Polyclonal Antibody

Catalog # **S19-63BR**

Lot # **Z2014-44**

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.