

Src monoclonal antibody (5A18)

Src is a member of the tyrosine kinase family that was originally identified as the oncogene of Rous sarcoma virus. The catalytic domain contains a tyrosine autophosphorylation site that increases kinase activity when phosphorylated. The Src family of tyrosine kinase is involved in a variety of receptor-induced signal transduction pathways such as immune recognition

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

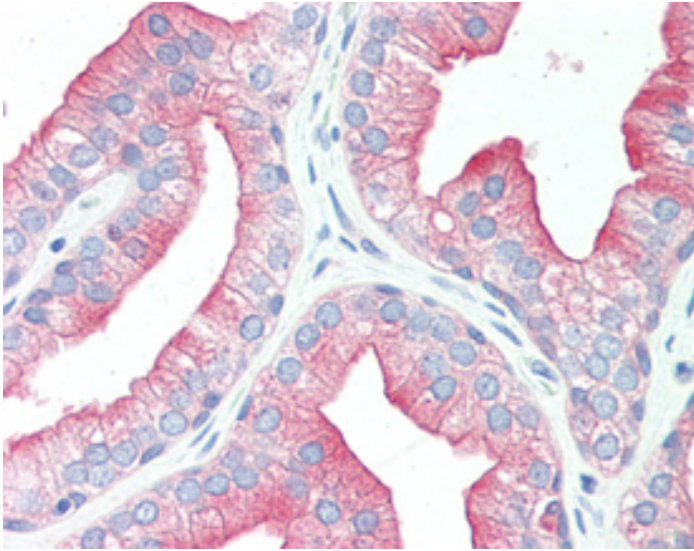
Ordering Information

[Order Online »](#)

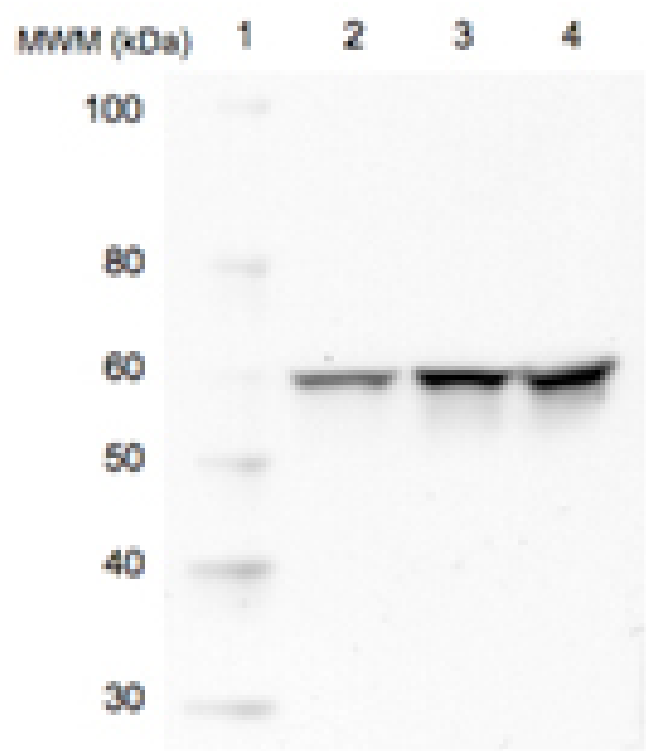
ADI-905-678-100	100µg
-----------------	-------

Manuals, SDS & CofA

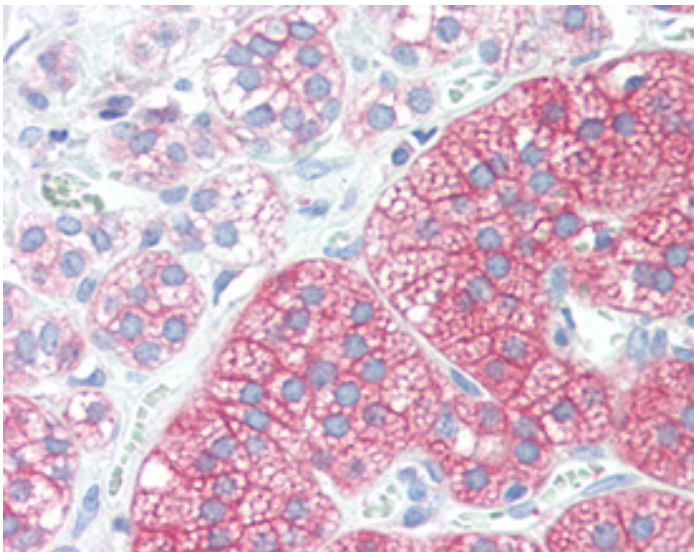
[View Online »](#)



IHC Analysis: Human prostate tissue stained with Src, mAb (5A18) at 20µg/ml.



Western blot analysis of Src using 20µg of whole cell lysate (Lane 2=HeLa, Lane 3=3T3, Lane 4=PC12) probed with with Src, mAb (5A18) at 1µg/ml.



IHC Analysis: Human adrenal tissue stained with Src, mAb (5A18) at 20µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proto-oncogene c-Src, pp60c-src

Application IHC (PS), WB

Application Notes Detects a band of ~60kDa by Western blot.

Clone 5A18

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.

Host Mouse

Immunogen Recombinant human full length c-Src.

Purity Detail Protein G affinity purified.

Recommendation
Dilutions/Conditions Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rat

UniProt ID P12931

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)