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 rec

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

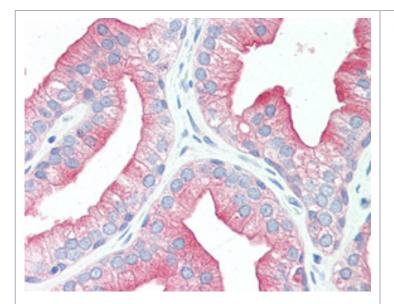
Order Online »

ADI-905-678-100

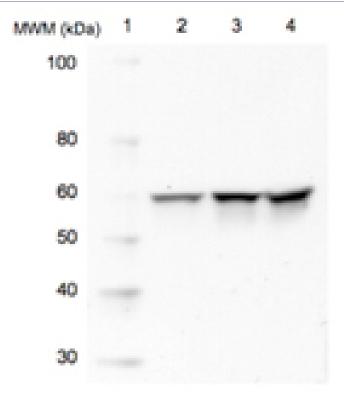
100µg

Manuals, SDS & CofA

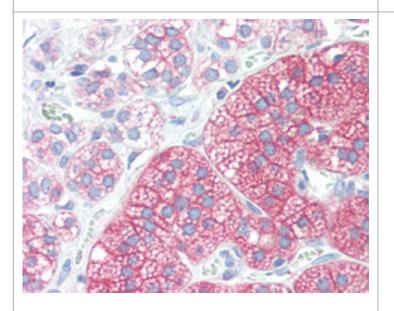
View Online »



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



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



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

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 NotesDetects 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

