

# Shc monoclonal antibody (11F6)

The Src homology 2 domain containing (Shc) adaptor protein links growth factor, cytokine, and antigen receptor stimulation to the Ras signaling pathway, and plays a critical role in insulin signaling, tumorigenesis, and hematopoietic cell growth and differentiation. Structurally, Shc is composed of an N-terminal phosphotyrosine binding (PTB) domain, a medial CH1 domain, and a C-terminal SH2 domain. Upon binding to activate cell-surface receptors, Shc is phosphorylated at three tyrosine residues (Tyr239, Tyr240, Tyr317) triggering recognition of Shc by SH2-containing proteins such as Grb2.

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

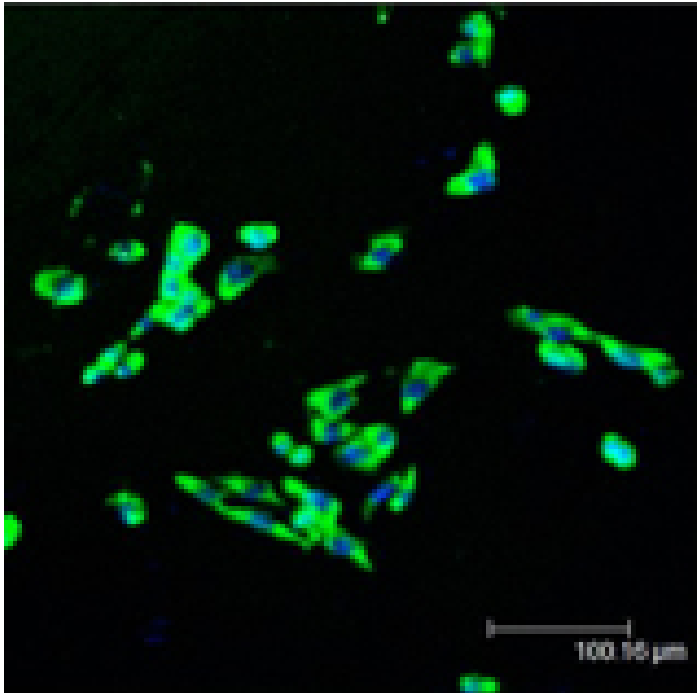
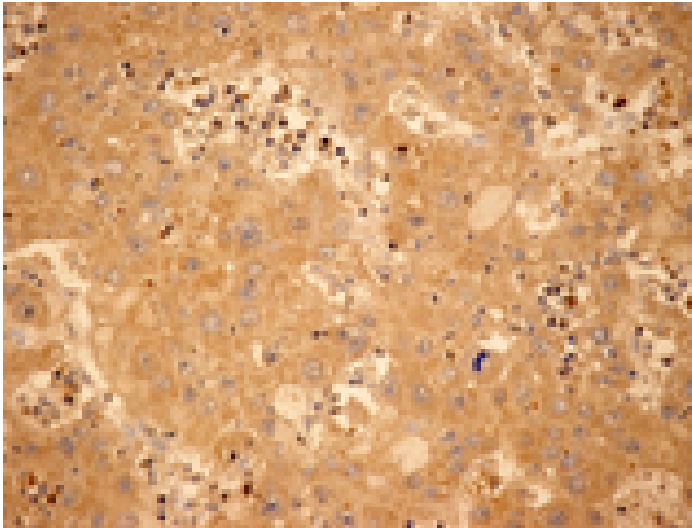
## Ordering Information

[Order Online »](#)

ADI-905-665-100	100µg
-----------------	-------

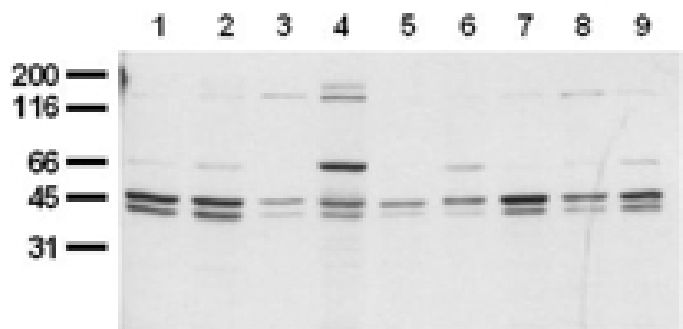
## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human hepatoma tissue with Shc mAb (11F6).

Immunofluorescent analysis (confocal) staining of V79 cells using Shc mAb (11F6) (green); nuclei are stained in blue pseudocolor using DRAQ5.



Western blot analysis of cancer cell lysates with Shc mAb (11F6); HeLa (1), HepG2 (2), HEK293 (3), SH-SY5Y (4), MDCK (5), PC12 (6), CMT93 (7), Neuro 2A (8), NIH-3T3 (9).

## Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Src homology 2 domain containing
Application	ELISA, ICC, IHC, WB
Application Notes	Detects bands of ~46kDa, ~52kDa and ~66kDa by Western blot.
Clone	11F6
Formulation	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of Shc.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Reconstitution	Reconstitute with 1ml water (15 minutes at room temperature).
Species Reactivity	Human, Mouse
UniProt ID	P29353



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

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

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

France  
Phone: +33 1 402 44 555  
[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)

This antibody is covered by our **Worry Free Guarantee**