

[pTyr⁶⁹⁵/Tyr⁶⁹⁹]Stat5 A/B monoclonal antibody (5G4)

Stats (Signal transducers and activators of transcription) are a family of cytoplasmic latent transcription factors that are activated to regulate gene expression in response to a large number of extracellular signaling polypeptides including cytokines, interferons, and growth factors. After phosphorylation by JAK tyrosine kinases, Stats enter the nucleus to regulate transcription of many different genes.

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

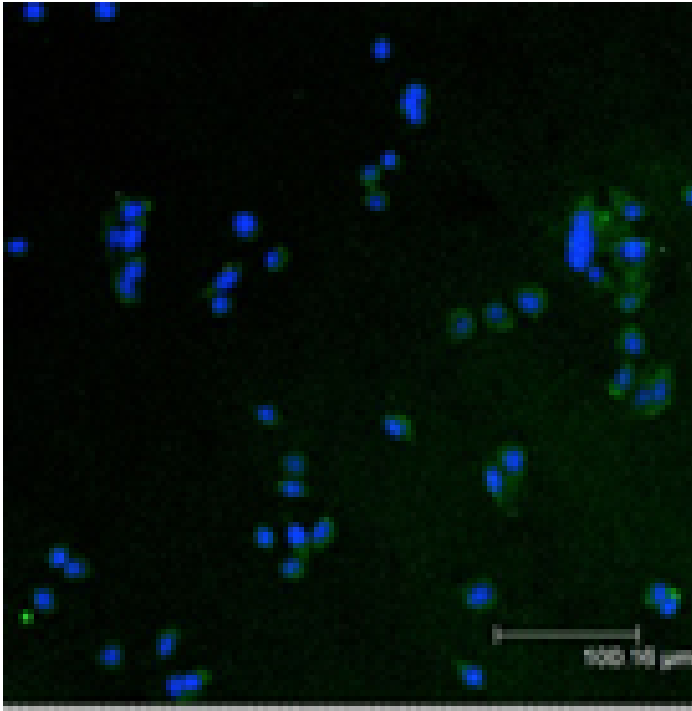
Ordering Information

[Order Online »](#)

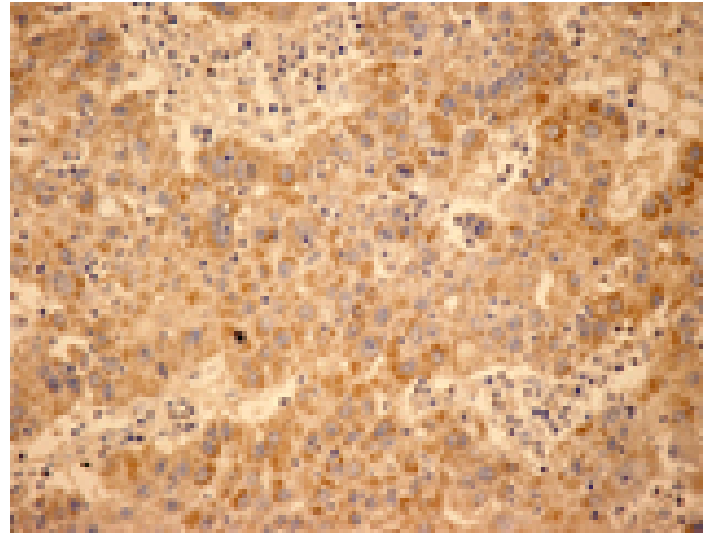
ADI-905-673-100	100µg
-----------------	-------

Manuals, SDS & CofA

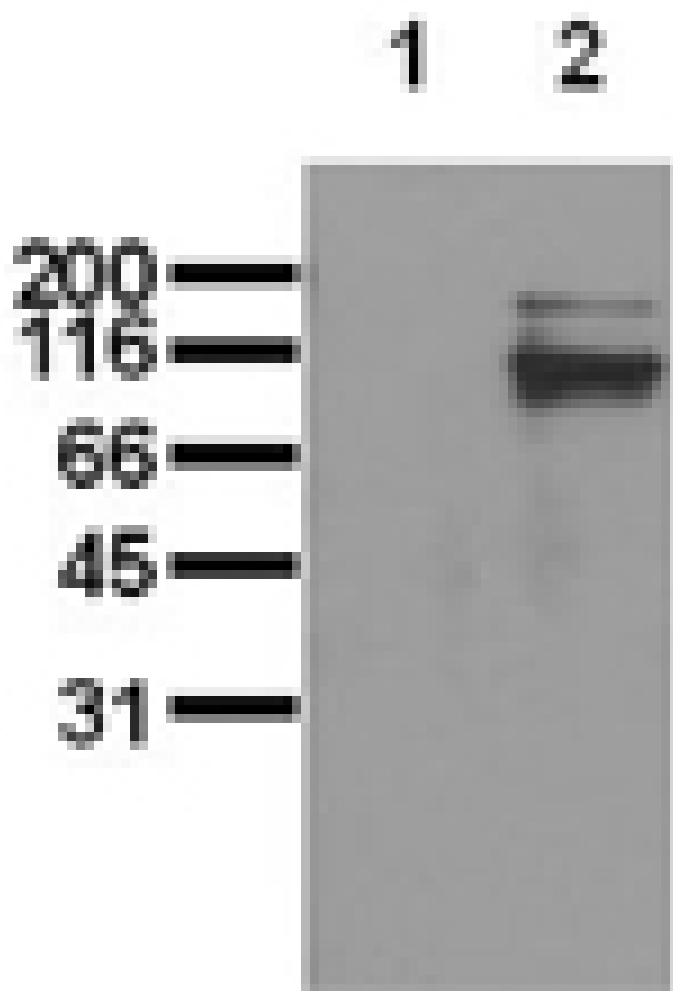
[View Online »](#)



Immunofluorescent analysis (confocal) staining of MCF-7 cells using Stat5A/B (pTyr694/699) mAb (5G4) (green); nuclei are stained in blue pseudocolor using DRAQ5.



Immunohistochemistry analysis of human hepatoma tissue with Stat5A/B (pTyr694/699) mAb (5G4).



Western blot analysis of untreated (1) and pervanadate treated (2) MCF-7 cells, probed with Stat5A/B (pTyr694/699) mAb (5G4).

Handling & Storage

Use/Stability Stable at -20°C up to 1 year.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Signal transducer and activator of transcription 5A/B

Application ELISA, ICC, IHC, WB

Clone 5G4

Formulation Liquid. In PBS containing 50% glycerol, PEG, sucrose, and 0.09% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of Stat5A/B phosphorylated at Tyr695/699.

Isotype IgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography purified.

Recommendation Dilutions/Conditions ELISA (0.05µg/ml)Immunohistochemistry (1:50)Western Blot (0.5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P42229 (STAT5A), P51692 (STAT5B)

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)