

# [pSer<sup>727</sup>]Stat3 monoclonal antibody (23G5)

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.

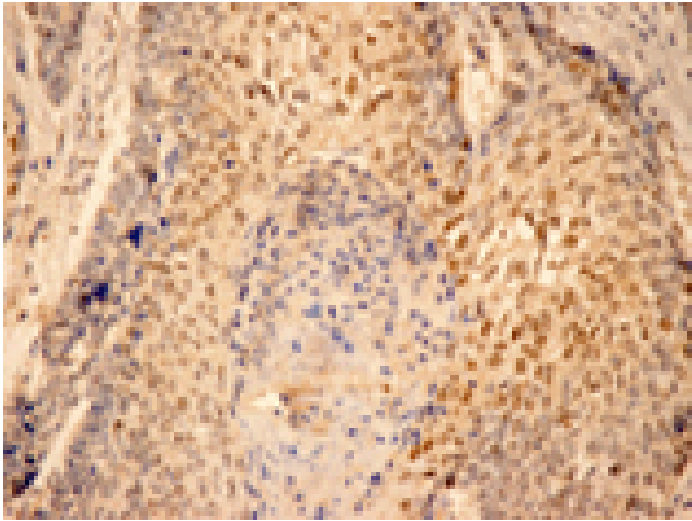
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

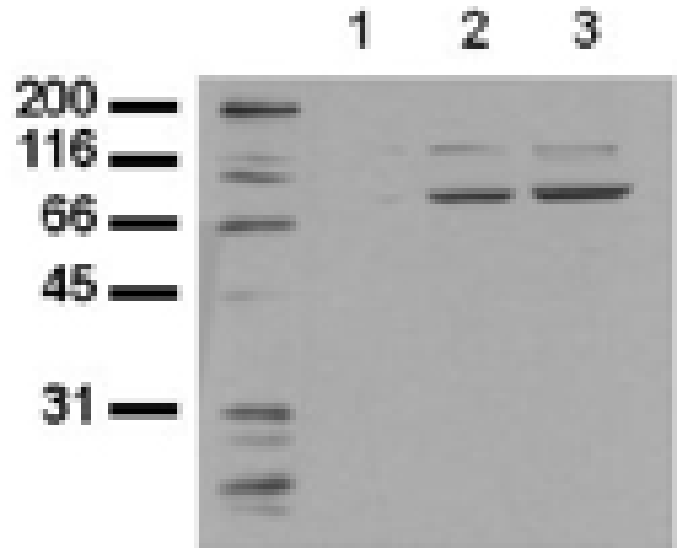
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human lung cancer tissue with Stat3 (pSer727) mAb, clone 23G5.



Western blot analysis of untreated (1), EGF stimulated (2), or pervanadate treated (3) A431 cell lysates with Stat3 (pSer727) mAb, clone 23G5.

## 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** Signal transducer and activator of transcription 3

**Application** ELISA, IHC, WB

**Application Notes** Detects a band of ~92kDa by Western blot.

**Clone** 23G5

**Formulation** Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to a portion of Stat3 phosphorylated at Ser727.

**Isotype** IgG1

**Purity Detail** Thiophilic adsorption and size exclusion chromatography purified.

**Recommendation Dilutions/Conditions** ELISA (0.1µg/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.

**Reconstitution** Reconstitute with 1ml water (15 minutes at room temperature).

**Species Reactivity** Dog, Human, Mouse, Rat

**UniProt ID** P40763

Last modified: May 29, 2024



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)