

# [pTyr<sup>705</sup>]STAT3 monoclonal antibody (9E12)

The STAT proteins serve as both cytoplasmic signal transducers and nuclear activators of transcription. STAT3 is activated by tyrosine phosphorylation at residue Y705 in cells treated with interleukin-6 or EGF. Activated STAT3 can bind to DNA either as homodimer or as heterodimer with STAT1.

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

## Ordering Information

[Order Online »](#)

ALX-804-180-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable at -80°C up to 1 year, at 4°C up to 3 months.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Signal transducer and activator of transcription 3
<b>Application</b>	ELISA, ICC, IHC, IP, WB
<b>Clone</b>	9E12
<b>Crossreactivity</b>	Does not cross-react with the non-phosphorylated form of STAT3 nor with unrelated tyrosine-phosphorylated proteins.
<b>Formulation</b>	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
<b>Host</b>	Mouse
<b>Immunogen</b>	Synthetic peptide corresponding to a portion of STAT3 phosphorylated at Tyr705 conjugated to KLH.
<b>Isotype</b>	IgG1
<b>Positive Control</b>	Included. (Prod. No. ALX-840-025)
<b>Purity Detail</b>	Thiophilic adsorption and size exclusion chromatography purified.

**Recommendation Dilutions/Conditions**

Western Blot (0.1µg/ml for HRPO/ECL detection; recommended blocking buffer CPPT: 10mM TRIS-HCl, pH 7.4, 0.5% (w/v) casein, 1% (w/v) PEG 4,000, 1% (w/v) polyvinylpyrrolidone, 0.1% (v/v) Tween 20, 150mM sodium chloride)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

**Specificity**

Recognizes the -AAPpYLKT-motif of activated STAT3 phosphorylated at Tyr<sup>705</sup>.

**Technical Info / Product Notes**

Replacement for **ADI-905-671**

**UniProt ID**

P40763

**Worry-free Guarantee**

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

**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)