

[pTyr⁷⁰⁵]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-804-1800	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
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, ICC, IHC, IP, WB
Clone	9E12
Crossreactivity	Does not cross-react with the non-phosphorylated form of STAT3 nor with unrelated tyrosine-phosphorylated proteins.
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 Tyr705 conjugated to KLH.
Isotype	IgG1
Positive Control	Included. (Prod. No. ALX-840-025)
Purity Detail	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 Tyr⁷⁰⁵.

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com