

Syk monoclonal antibody (SYK-01)

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

Citations: 8

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ICC, IHC (PS), IP, WB
Clone	SYK-01
Formulation	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
Host	Mouse
Immunogen	Recombinant fragment (aa 5-360) of human Syk.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 5µg/ml)Western Blot (1-2µg/ml, 60min.)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P43405
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