

Phosphotyrosine monoclonal antibody (PY20) (agarose immobilized)

Covalently coupled to 4% cross-linked agarose beads. This product is optimized for binding of phosphotyrosine containing proteins with minimal non-specific binding.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Immunoaffinity chromatography, IP
Clone	PY20
Crossreactivity	No cross-reactivity with phosphoserine or phosphothreonine.
Formulation	50% slurry in PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Phosphotyrosine.
Isotype	IgG2b
Source	Purified from mouse ascites.
Species Reactivity	Species independent
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