

# Phosphotyrosine monoclonal antibody (PY20)

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SA240-1000	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	ICC, IHC, IP, WB
<b>Clone</b>	PY20
<b>Crossreactivity</b>	Does not cross-react with phosphoserine or phosphothreonine.
<b>Formulation</b>	Liquid. In PBS containing 50% glycerol and 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Phosphotyrosine.
<b>Isotype</b>	IgG2b
<b>Purity Detail</b>	Purified.
<b>Source</b>	Purified from mouse ascites.
<b>Species Reactivity</b>	Species independent
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@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)