

# HSP90 monoclonal antibody (3G3)

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

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-079-R400	400µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Before use, centrifuge suspension (4'000 x g for 15 min.) and discard supernatant. Resuspend pellet in 200µl PBS, pH 7.2. If desired, add 0.05% sodium azide. Stable for 6 months after PBS resuspension. Removal of residual ammonium sulfate is not usually necessary for IP procedure; for other uses it may be removed by dialysis or equivalent method.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Heat shock protein 90
<b>Application</b>	IF, IHC, IP, WB
<b>Clone</b>	3G3
<b>Formulation</b>	Liquid. Contains no preservatives.
<b>Host</b>	Mouse
<b>Immunogen</b>	Purified mouse HSP90 (heat shock protein 90).
<b>Isotype</b>	IgM
<b>Purity Detail</b>	Ammonium sulfate/ascites suspension.
<b>Recommendation Dilutions/Conditions</b>	Immunofluorescence (1:200)Immunohistochemistry (1:200)Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Chicken, Human, Mouse, Rainbow trout, Rat
<b>UniProt ID</b>	P07901 (HSP90α), P11499 (HSP90β)

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)