

HSP90 monoclonal antibody (3G3)

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

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Heat shock protein 90
Application	IF, IHC, IP, WB
Clone	3G3
Formulation	Liquid. Contains no preservatives.
Host	Mouse
Immunogen	Purified mouse HSP90 (heat shock protein 90).
Isotype	IgM
Purity Detail	Ammonium sulfate/ascites suspension.
Recommendation Dilutions/Conditions	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.
Species Reactivity	Chicken, Human, Mouse, Rainbow trout, Rat
UniProt ID	P07901 (HSP90α), P11499 (HSP90β)

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)