

# Metallothionein (human) clonal antibody (N11-G)

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 12 months after receipt when stored as recommended.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	ELISA, WB
<b>Clone</b>	N11-G
<b>Formulation</b>	Liquid. In 20mM TRIS/HCl, pH 8.0, containing 10mg/ml BSA and 0.05% sodium azide.
<b>Host</b>	Rabbit
<b>Immunogen</b>	Synthetic peptide corresponding to the N-terminal sequence of human metallothionein (covering isoforms 1A, B, E, F, G, H, I, L, M, Q, S, Y, 2A).
<b>Isotype</b>	IgG
<b>Purity Detail</b>	Major clone of IgGs obtained from immunoaffinity purified immunoglobulins corresponding to immunogenic peptide.
<b>Recommendation Dilutions/Conditions</b>	ELISA (1:5,000-1:10,000)Western Blot (1:1,000; Wash buffer: 1x TRIS buffered saline (TBS) containing 0.1% Triton X-100, Blocking buffer: 1x TBS containing 0.1% Triton X-100 and 8% nonfat dry milk. Incubate the membrane with antibody diluted in blocking buffer for 2 hours at room temperature)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human

**Clonal Rabbit Antibody** is a pure homogenous fraction of rabbit immunoglobulin (IgG) which corresponds to a unique linear epitope on the target protein. This epitope is designed after a detailed analysis of the antigen structure, regulatory post-translational modifications, protein/protein interaction domain and protein folding. The chosen sequence (peptide) is synthesized and the rabbit is immunized. The clone is then separated from the crude antiserum by “*in vitro* cloning technology”. The resulting product is monospecific and characterized by superior specificity, affinity and avidity.

**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)