

Metallothionein monoclonal antibody (UC1MT)

Metallothioneins (MTs) are a class of ubiquitous, low molecular weight, cysteine-rich, metal binding proteins that function in the metabolism of zinc and copper. MT protects against exposure to toxic metals such as mercury and cadmium as well as a variety of other stress conditions. There are four mammalian MT isoforms: MT-I, -II, -III, -IV, each of which is encoded by individual genes. MT expression is induced by heavy metal cations, starvation, elevated temperature, lipopolysaccharide, inflammatory cytokines, free radicals, and ionizing radiation.

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

Citations: 12

[View Online »](#)

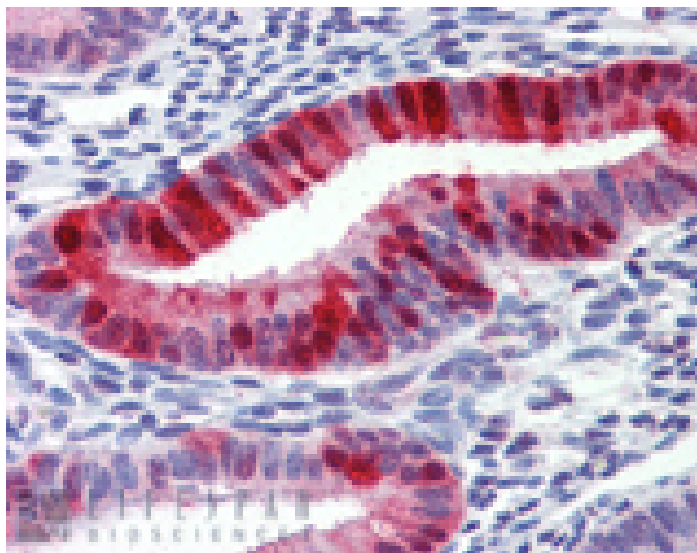
Ordering Information

[Order Online »](#)

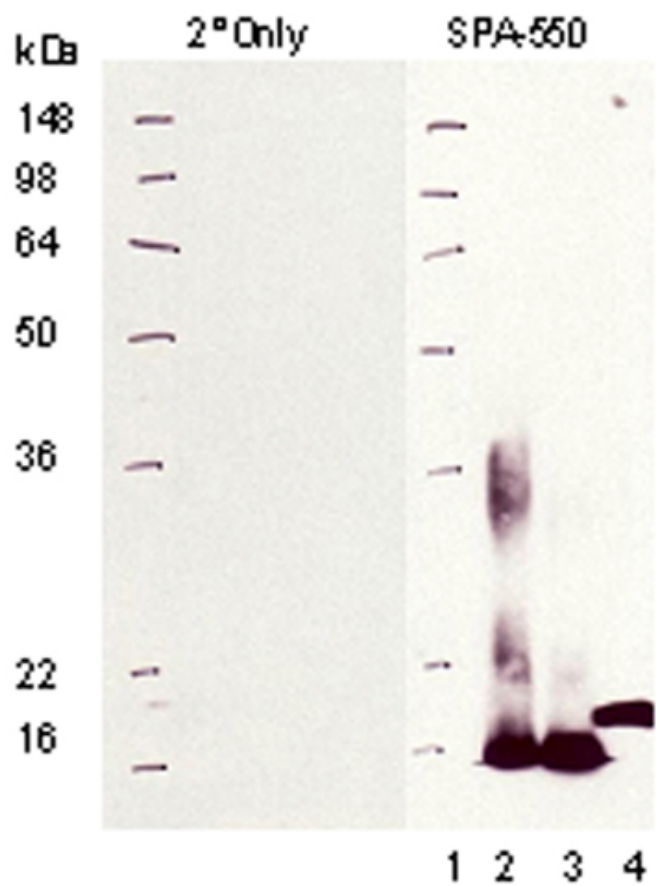
ADI-SPA-550-D	50µg
ADI-SPA-550-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human uterus tissue stained with Metallothionein, mAb (UC1MT) at 10µg/ml.



Western blot analysis of Prod. No. ADI-SPA-550: Lane 1: MW marker, Lane 2: MT I, Lane 3: MT II, Lane 4: Mummichug CdCl₂.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MT1A, Metallothionein MGC32848, MT1S, MTC, metallothionein 1A (functional)
Application	ELISA, IHC (PS), WB
Application Notes	Detects a band of ~15kDa by Western blot.
Clone	UC1MT
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Cross-linked rabbit liver Metallothionein I and II.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Fish, Horse, Human, Mouse, Rabbit, Rat

Worry-free Guarantee

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



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)

Last modified: May 29, 2024