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: 11

View Online »

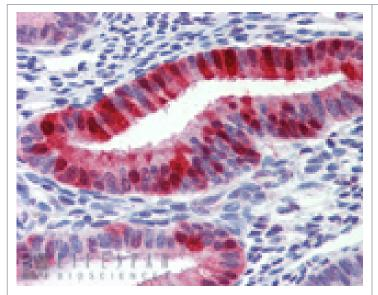
Ordering Information

Order Online »

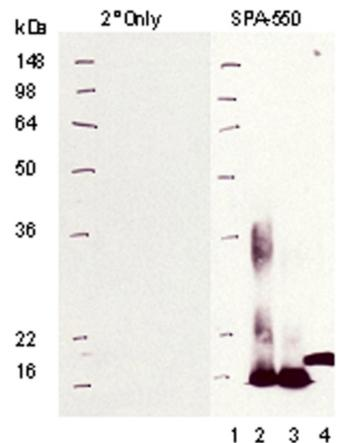
ADI-SPA-550-D	50µg
ADI-SPA-550-F	200μg

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.

IgG1 Isotype

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.

Purified from ascites. Source

Species Reactivity Fish, Horse, Human, Mouse, Rabbit, Rat

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

ENZO LIFE SCIENCES, Phone: 800.942.0430 info-Last modified: May 29, 2024 European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com