

# Cytochrome c monoclonal antibody (7H8.2C12)

Cytochrome c is an electron transport protein from mitochondria. A primary means of caspase activation by p53 is through induction of cytochrome c release from mitochondria. Cleavage of Bid, a Bcl 2 family member, by caspase-8 produces a truncated tBid molecule which triggers mitochondria release of cytochrome c which binds to the adaptor protein Apaf-1 and procaspase-9 to form a cytosolic apoptosome complex. This event results in activation of procaspase-9 and its subsequent cleavage and activation of caspase-3 leading to the degradation of a series of cellular substrates and subsequently cell death.

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

Citations: 1

[View Online »](#)

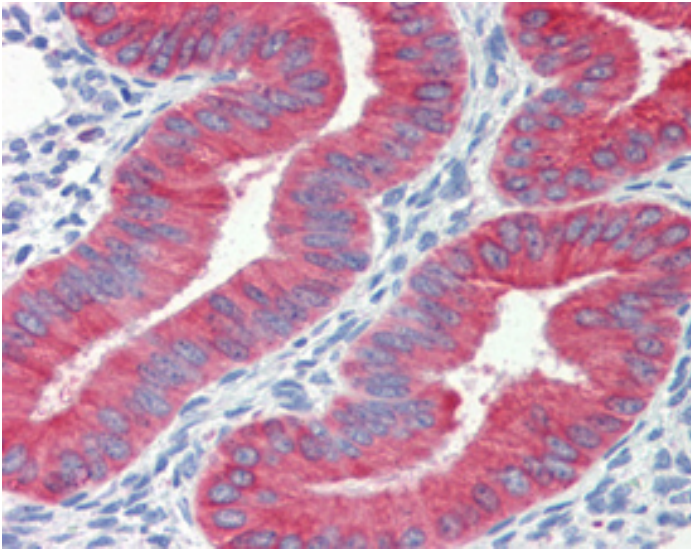
## Ordering Information

[Order Online »](#)

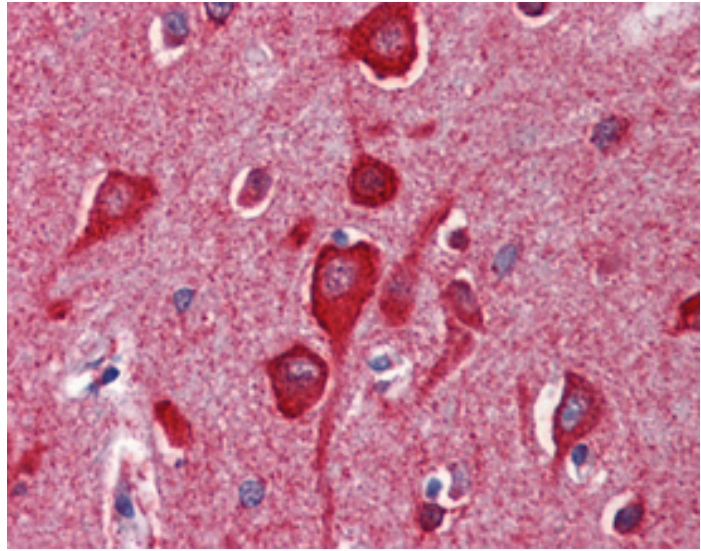
ADI-AAM-175-E	100µg
---------------	-------

Manuals, SDS & CofA

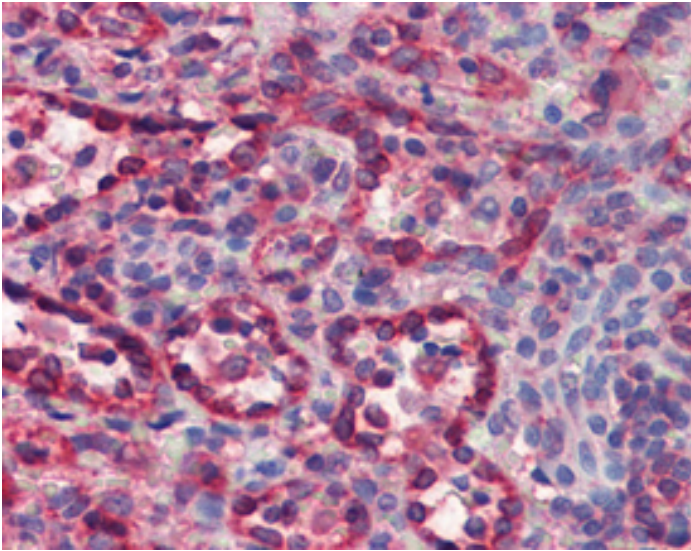
[View Online »](#)



IHC Analysis: Human uterus tissue stained with Cytochrome c, mAb (7H8.2C12) at 5µg/ml.



IHC Analysis: Human brain, cortex tissue stained with Cytochrome c, mAb (7H8.2C12) at 5µg/ml.



IHC Analysis: Human spleen tissue stained with Cytochrome c, mAb (7H8.2C12) at 5µg/ml.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cyt c

**Application** IHC (PS), WB

**Application Notes** Detects a band of ~14kDa by Western blot.

**Clone** 7H8.2C12

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**Host** Mouse

**Immunogen** Horse Cytochrome c .

**Isotype** IgG2b

**Purity Detail** Protein G affinity purified.

**Recommendation** Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available  
**Dilutions/Conditions** for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from mouse ascites.

**Species Reactivity** Bovine, Guinea pig, Hamster, Human, Monkey, Mouse, Rabbit, Rat, Sheep

**UniProt ID** P00004



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 42 44 55 55  
[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)