

Cytochrome c monoclonal antibody (6H2.B4)

Cytochrome c, which is a critical component of the electron transport chain, is a component of the functional apoptosome. Upon release from the mitochondria, cytochrome c binds to the adaptor proteins 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](#).

Manuals, SDS & CofA

[View Online »](#)

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 ELISA, IP

Application Notes Estimated molecular weight of ~13kDa.

Clone 6H2.B4

Formulation Liquid. In PBS containing 0.1% sodium azide.
Concentration 1mg/ml.

Host Mouse

Immunogen Native rat Cytochrome c.

Isotype IgG1

Purity Detail Protein G affinity purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rat

UniProt ID P62898

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

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

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

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

Last modified: May 29, 2024