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.

Ordering Information

Order Online »

ADI-905-634-100 100μg

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.

lgG1

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655
infofr@enzolifesciences.com

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

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