Cyclophilin A polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 32

View Online »

Ordering Information

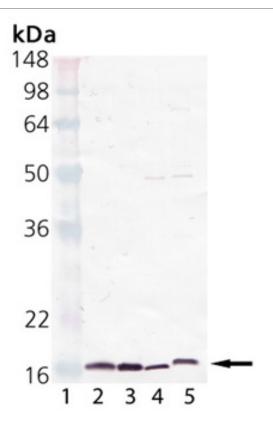
Order Online »

BML-SA296-0100

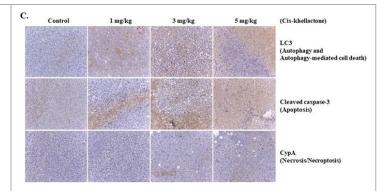
100µl

Manuals, SDS & CofA

View Online »

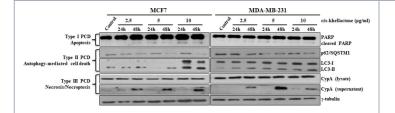


Western Blot analysis of Cyclophilin A, pAb . Lane 1: M.W. Marker, Lane 2: BML-SE105, Lane 3: Jurkat Cell Lysate, Lane 4: Rat Brain Extract, Lane 5: Mouse Brain Extract



In vivo assessment of cis-khellactone anti-tumor activity in a murine model(A) Change in tumor volume after ciskhellactone treatment of a nude mouse. Tumor-bearing mice were injected with MDA-MB-231 cancer cells as mentioned in Materials and Methods. When tumors were approximately 50 to 100 mm3 in volume, mice in each treatment group were intravenously injected via the tail vein with cis-khellactone (at a dose of 1, 3, or 5 mg/kg) once every 3 days. Control groups received only normal saline (*P<0.01, n = 5, Student's t test). (B) Representative images of H&E staining of five major organs (heart, lung, liver, spleen, and kidney) and tumor. Mice were sacrificed at 30 days after the initial drug administration and tissue samples were immediately collected. Scale bar = 140 µm. (C) Induction of three PCD in xenograft mice tumor model with MDA-MB-231 cells were tested by employing Immunohistochemistry (IHC). Three PCD were checked with cleaved caspase 3 for apoptosis, LC3 for autophagic cell death, and CypA for necrosis/necroptosis.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Anti-cancerous effect of cis-khellactone from Angelica amurensis through the induction of three programmed cell deaths. *Oncotarget* (2018)



Effect of cis-khellactone on three types of PCD (apoptosis, autophagy-mediated cell death, and necrosis/necroptosis) in MCF7 and MDA-MB-231 cancer cellsMCF7 and MDA-MB-231 cells were cultured and treated with 2.5, 5, and 10 μg/ml for 24 or 48 h, while control cells were treated with DMSO alone. Cell lysates were examined by Western blot analysis by using following corresponding biomarkers or regulatory proteins of the different types of cell death; PARP for apoptosis, p62 and LC3 for autophagy, and CypA for necrosis/necroptosis; γ-tubulin was used as an internal control. The result of Western blot was quantified by using ImageJ program.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Anti-cancerous effect of cis-khellactone from Angelica amurensis through the induction of three programmed cell deaths.

Oncotarget (2018)

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CyPA, Peptidyl-prolyl cis-trans isomerase A

Application ICC, WB

Formulation Liquid. Neat serum.

Host Rabbit

Immunogen Recombinant human cyclophilin A conjugated to KLH.

Source Whole rabbit serum.

Species Reactivity Human, Mouse, Rat

UniProt ID P62937

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

.

Last modified: May 29, 2024

