

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

Anti-Caspase-9 Monoclonal Antibody (clone LAP6-96)

Catalog #: 2308-MC-050

Size: 50 µg

Description: Caspases are key effectors of programmed cell death. They are synthesized as inactive proenzymes which are activated by cleavage at a specific aspartate residue to form two subunits. These subunits are normally linked together by a linker which may be involved in the regulation of the different caspases. Caspase-9 is a member of the CED-3 family and bear high similarity to caspase-3. Procaspase-9 can be activated by either caspase-3 or granzyme B, although they cleave the proenzyme to subunits of different sizes. Cleavage by granzyme B produces an active enzyme which is capable of cleaving PARP. Also, the ability of caspase-3 to activate caspase-9 seems to suggest that caspase-9 is further downstream of caspase-3 and may be involved in later changes in cells observed undergoing apoptosis.

Physical State: This mouse monoclonal antibody is provided as purified IgG in phosphate buffer saline with 0.08% sodium azide. The final antibody concentration is 1 mg/ml.

Immunogen: Recombinant human caspase-9 protein.

Isotype: IgG₁

Specificity: The antibody detects human caspase-9.

Storage: Store at -20°C in working aliquots to avoid freeze-thaw cycles.

Applications: For western blotting, a starting concentration of 1-5 µg/ml is suggested. Empirical determination will be required for optimal results.



Figure 1. Western blot analysis using Caspase-9 antibody on MCF-7 cells negative (-) and positive (+) for caspase-3 and showing the proenzyme form of caspase-9 and one of the cleavage products after treatment with thapsigargin for 48 hours.

© 2001 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

v20716

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com

References:

Cohen, G.M. 1997. Caspases: the executioners of apoptosis. *Biochem J.* **326**:1-16.

Stennicke, H.R., Q.L. Deveraux, E.W. Humke, J.C. Reed, V.M. Dixit, and G.S. Salvesen 1999. Caspase-9 can be activated without proteolytic processing. *J. Biol. Chem.* **274**:8359-8362.

Kuida, K. 2000. Caspase-9. *Int. J. Biochem. Cell Biol.* **32**:121-124.

**Anti-Caspase-9
Monoclonal Antibody**

Catalog #: 2308-MC-100

Storage: -20°C

TREVIGEN[®]

1-800-873-8443