

## Caspase-8 Monoclonal Antibody [3B10]

(Cat. No. A-0566)

### Background

Caspase-8 and the sequential activation of caspases such as caspase-1, 3, 6, and 7, is important to the processes of cell apoptosis, or programmed cell death. The caspase-8 protein is thought to interact with Fas-interaction protein FADD and may also play a role in neurodegenerative diseases such as Huntington disease.

### Description

Rat monoclonal antibody to Caspase-8, Clone 1G12

### Formulation

Liquid. In PBS containing 0.02% sodium azide.

### Immunogen

Recombinant mouse caspase-8 p18 subunit

### Specificity

Mouse

### Isotype

IgG1

### Purification

Protein G- affinity purified.

### Storage

Store at +4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

### Application

ELISA, FC, ICC, WB

### Ordering Information

#### Products

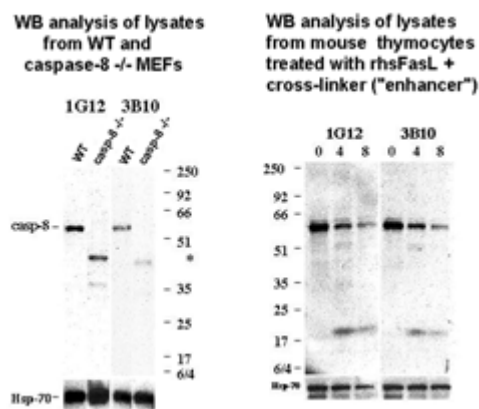
Caspase-8 Monoclonal Antibody [3B10]

#### Size

100 µg

#### Cat. No.

A-0566-100



**Fig. 1.** Caspase-8 Monoclonal Antibody [3B10] and [1G12] both detect pro-caspase-8 in MEFs from WT mice, but not in MEFs from caspase-8<sup>-/-</sup> mice. Several smaller bands detected in the caspase-8<sup>-/-</sup> MEFs, correspond to truncated forms of caspase-8 made in the caspase-8<sup>-/-</sup> mice since only exons 1 and 2 of mouse caspase-8 were deleted in these knock-out mice and not the region encoding the p18 subunit.