

# Granzyme B Monoclonal Antibody [B18.1], FITC Conjugated

(Cat. No. A-0526)

## Background

Granzymes B is serine proteases family member that is found in the granules of cytotoxic T cells (CTLs) and natural killer (NK) cells. Once a target cell is recognized, CTLs and NK cells secrete GrB and other granule proteins, including perforin, into the immunological synapse. GrB cleaves and activates, or inactivates, various protein substrates, resulting in the eventual apoptotic demise of the target cell. GrB apoptosis is also activated by adenovirus, which can effectively replace perforin. Extracellular granzyme B has been associated in the pathogenesis of vascular leakage, atherosclerosis aneurysm, chronic wound healing, and skin aging.

## Description

Mouse monoclonal antibody to Granzyme B, clone B18.1 (FITC conjugated)

## Formulation

Liquid. In PBS containing 0.02% sodium azide.

## Immunogen

Recombinant human granzyme B (aa 7-227).

## Specificity

Human

## Isotype

IgG1

## Purification

Serum

## Storage

Store at 4°C. Avoid multiple freeze/thaw cycles.

## Application

1µl of a 1mg/ml stock solution were used to stain 200'000 cells in a sample volume of 50µl, FC, ICC, IHC



**Fig. 1.** Detection of granzyme B in permeabilized human CD8+ T lymphocytes (filled profile). Isotype control in combination with a secondary FITC-labeled antibody is shown as control.

## Ordering Information

Product	Size	Cat. No.
Granzyme B Monoclonal Antibody [B18.1], FITC Conjugated	100 µl	A-0526-100

*This product is for research purposes only. Not intended for use in diagnostic procedures.*

Epigentek Group Inc. • USA Tel: 1-877-374-4368 • Int. Tel: 1-718-484-3990 • Fax: 1-718-484-3990 • [info@epigentek.com](mailto:info@epigentek.com) • [www.epigentek.com](http://www.epigentek.com)

2015-10-27-A-0526