

Perforin Monoclonal Antibody [CB5.4]

(Cat. No. A-0525)

Description

Rat monoclonal antibody to Perforin, Clone CB5.4

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant mouse perforin (aa 98-534)

Specificity

Mouse

Isotype

IgG2a

Purification

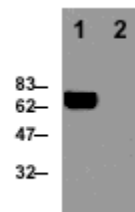
Serum

Storage

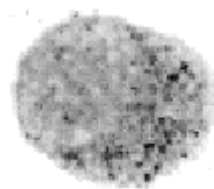
Store at +4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

Application

ICC, IHC, IP, WB



▲ **Fig. 1.** Detection of perforin in the mouse T cell clone B6.1. (lane 1). Absence of perforin in the mouse fibroblast cell line NIH.3T3 (lane 2). The protein migrates as a 66 kDa species.



▲ **Fig. 2.** Detection of perforin in an individual mouse T cell. Perforin is located in cytoplasmic granules.

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