## Bak (human) monoclonal antibody (TC100)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

**Ordering Information** 

Order Online »

BML-SA298-0050

50µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name BCL2L7

**Application** Flow Cytometry, ICC, WB

Clone TC100

**Formulation** Liquid. In PBS containing sodium azide and gelatin.

**Host** Mouse

Immunogen Recombinant N-terminal human bak.

lsotype lgG2a

Purity Detail Purified.

Species Reactivity Human

UniProt ID Q16611

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

.

Last modified: May 29, 2024

