

# 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.                                      |
| Isotype              | IgG2a                                                                  |
| Purity Detail        | Purified.                                                              |
| Species Reactivity   | Human                                                                  |
| UniProt ID           | Q16611                                                                 |
| Worry-free Guarantee | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> . |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)