

# TCR C $\beta$ 1 (human), mAb (JOVI-1)

## Ordering Information

[Order Online »](#)

ANC-101-040	FITC - 120 tests
-------------	------------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Flow Cytometry

Application Notes Functional Application (activates T cells expressing TCR Cβ1).

Clone JOVI-1

Host Mouse

Immunogen Human HA1.7 TCRβ chain expressed on transgenic mouse cells.

Isotype IgG2a

Species Reactivity Human

**Technical Info / Product Notes** For the Original Manufacturer's Data Sheet please click on the corresponding product number:

[ANC-101-020](#)

[ANC-101-030](#)

[ANC-101-040](#)

[ANC-101-050](#)

[ANC-101-060](#)

[ANC-101-820](#)



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)