

# BEST1 Antibody

## Ordering Information

[Order Online »](#)

PSC-15-355-R100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** IF, WB

**Formulation** Liquid. PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

**Host** Rabbit

**Immunogen** Recombinant fusion protein containing a sequence corresponding to amino acids 376-585 of human BEST1 (NP\_004174.1).

**Isotype** IgG

**Purity Detail** Affinity purification

**Species Reactivity** Human, Mouse

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).

**UniProt ID** O76090



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)