

# Oct4 monoclonal antibody (03)

This monoclonal antibody specifically detects Oct4, a transcription factor essential for maintaining the pluripotency and self-renewal of embryonic stem cells. It is widely used in stem cell research and developmental biology.

Oct4 (Octamer-binding transcription factor 4) is a POU-domain transcription factor crucial for maintaining the undifferentiated state of embryonic stem cells. It is expressed in the inner cell mass of blastocysts and in induced pluripotent stem cells (iPSCs). Dysregulation of Oct4 has been implicated in various cancers and is used as a marker for pluripotency.

This antibody is covered by our [Worry-Free Guarantee](#).

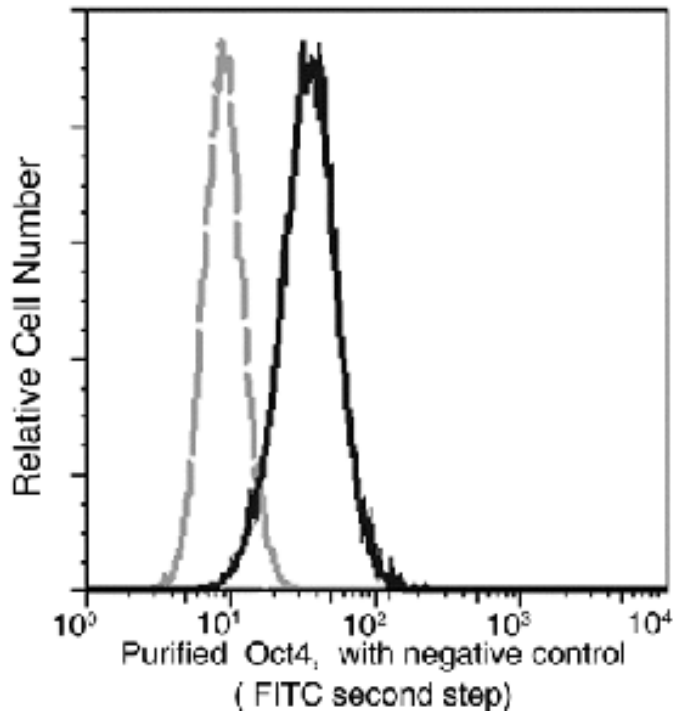
## Ordering Information

[Order Online »](#)

ENZ-ABS1063-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Oct4 in human embryonic stem (ES) cells. Human embryonic stem (ES) cells were fixed and permeabilized with BD Cytfix/Cytoperm™ Fixation/Permeabilization buffer (BD Pharmingen™ Cat. No. 554714), and stained with Purified Mouse anti-Oct4 monoclonal antibody, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

## Handling & Storage

**Handling** Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Octamer-binding protein 4, POU domain class 5 transcription factor 1, POU5F1, Oct-3, Octamer-binding transcription factor 3, OTF-3

**Application** Flow Cytometry

**Clone** 03

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to the C-terminus of human Oct4.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation** Flow Cytometry (1:25-1:100)  
**Dilutions/Conditions** Suggested dilutions/conditions may not be available for all applications.  
Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



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-be@enzolifesciences.com)

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