# 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**

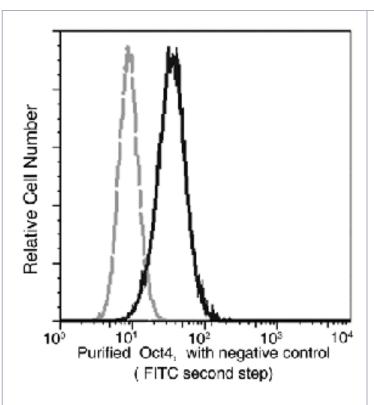
Order Online »

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 Cytofix/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 Dilutions/Conditions** Flow Cytometry (1:25-1:100)

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

•

Last modified: March 28, 2025



info-

eu@enzolifesciences.com