

# Aldolase C

## monoclonal antibody

### (AT2E11)

Aldolase C is a member of the class I fructose-biphosphate aldolase family. This protein is a glycolytic enzyme that catalyzes the reversible aldol cleavage of fructose-1, 6-biphosphate and fructose 1-phosphate to dihydroxyacetone phosphate and either glyceraldehyde-3-phosphate or glyceraldehydes respectively. It is expressed specifically in the hippocampus and Purkinje cells of the brain.

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

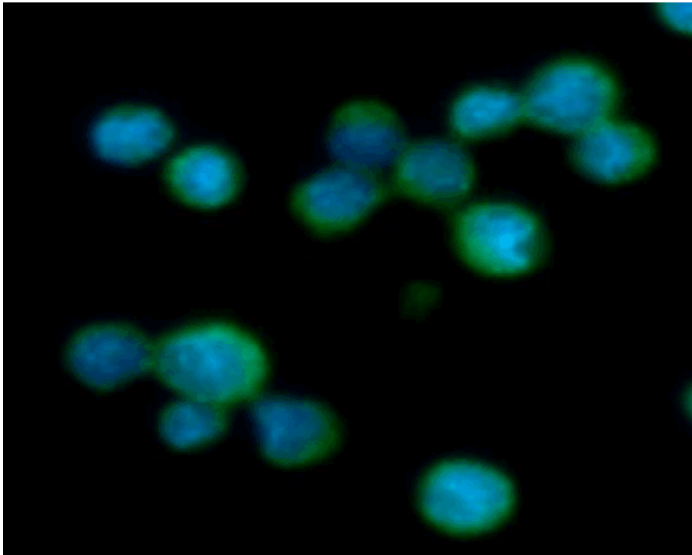
#### Ordering Information

[Order Online »](#)

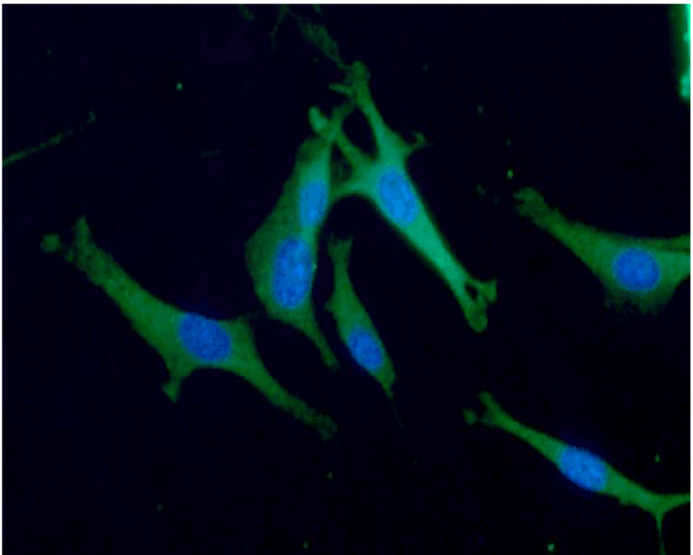
ENZ-ABS761-0100	100µl
-----------------	-------

#### Manuals, SDS & CofA

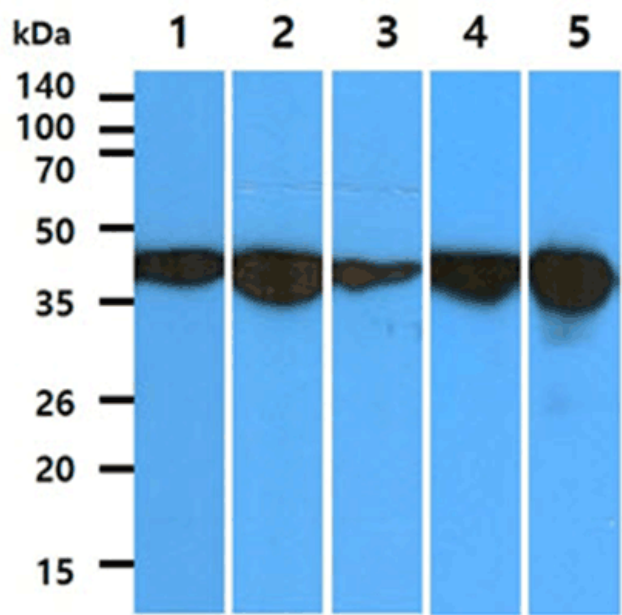
[View Online »](#)



IF analysis of ALDOC in Jurkat cell line, stained with DAPI (Blue) for nucleus staining and monoclonal anti-human ALDOC antibody (1:100) with goat anti-mouse IgG-Alexa fluor 488 conjugate (Green).



IF analysis of ALDOC in U87MG cell line, stained with DAPI (Blue) for nucleus staining and monoclonal anti-human ALDOC antibody (1:100) with goat anti-mouse IgG-Alexa fluor 488 conjugate (Green).



The cell line and tissue lysates (40µg) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with the ALDOC antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system. Lane 1: PC3 cell lysate, Lane 2: U87MG cell lysate, Lane 3: WiDr cell lysate, Lane 4: Jurkat cell lysate, Lane 5: Mouse brain tissue lysate.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	ALDC, Fructose biphosphate aldolase C, Brain-type aldolase
<b>Application</b>	ELISA, ICC, IF, WB
<b>Clone</b>	AT2E11
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human Aldolase C (aa 1-364) purified from E. coli.
<b>Isotype</b>	IgG2ak
<b>Purity Detail</b>	Protein A affinity purified
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	P09972
<b>Worry-free Guarantee</b>	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-fr@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)