

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](#).

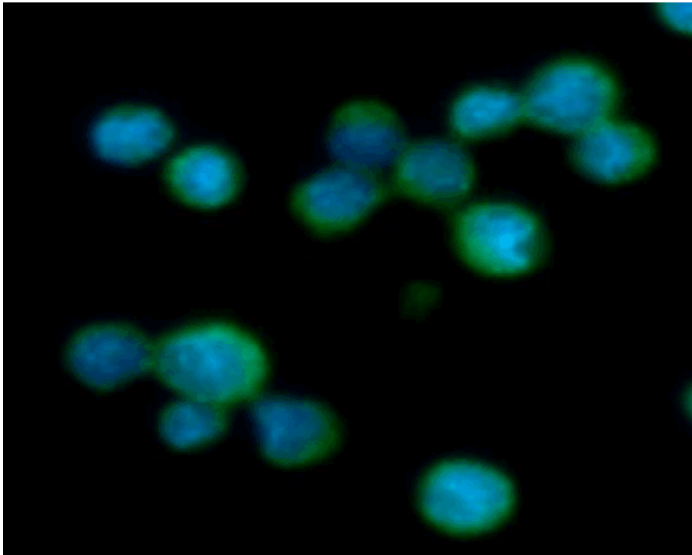
Ordering Information

[Order Online »](#)

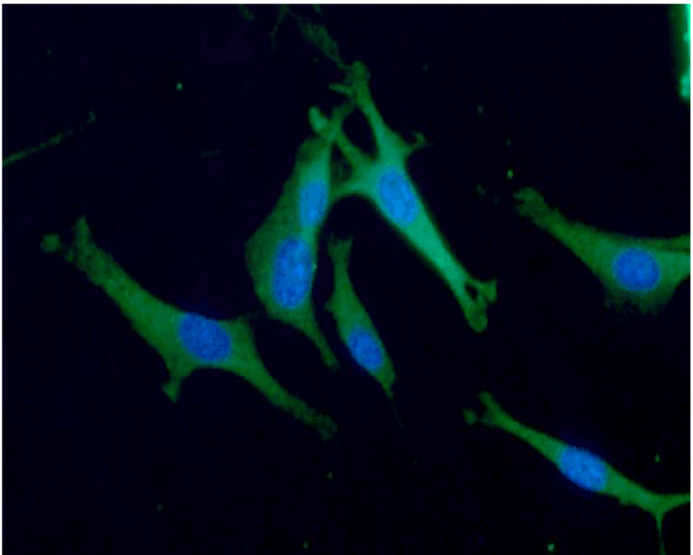
| | |
|-----------------|-------|
| ENZ-ABS761-0100 | 100µl |
|-----------------|-------|

Manuals, SDS & CofA

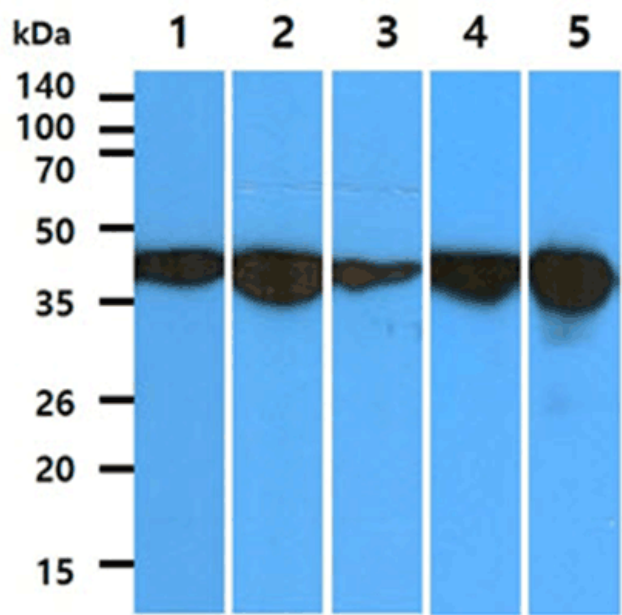
[View Online »](#)



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

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