

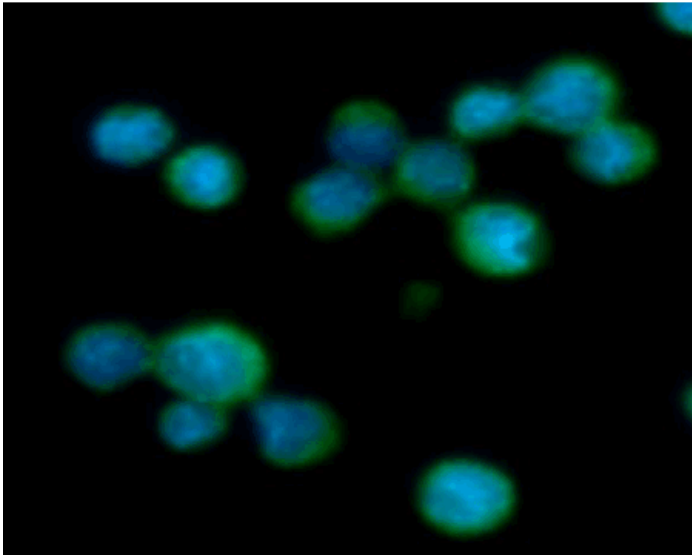
Aldolase C monoclonal antibody (AT2E11)

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.

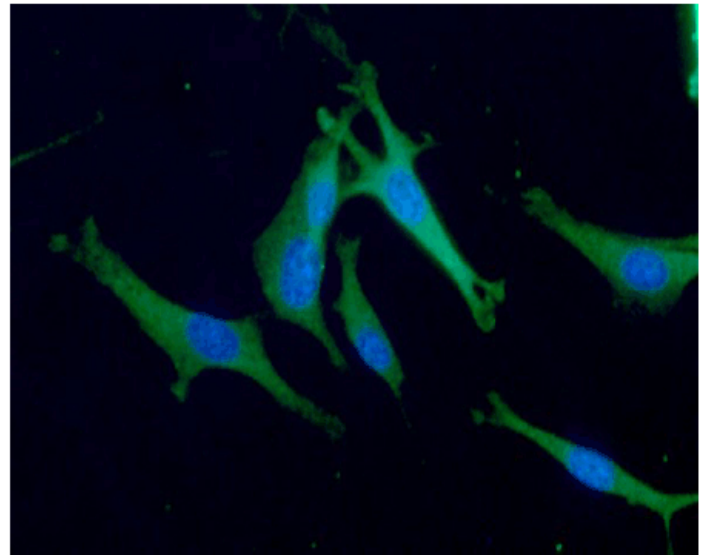
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

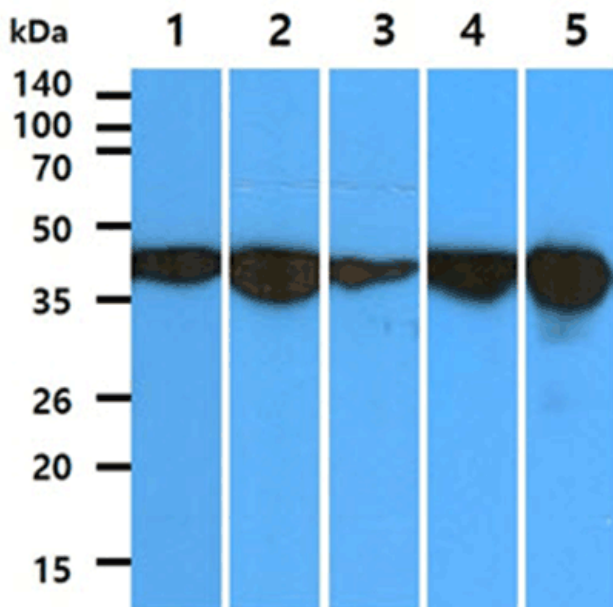
[View Online »](#)



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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com