

DLK1 (mouse) monoclonal antibody (PF183E)

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

Ordering Information

[Order Online »](#)

ALX-804-7430	50µg
--------------	------

Manuals, SDS & CofA

[View Online »](#)

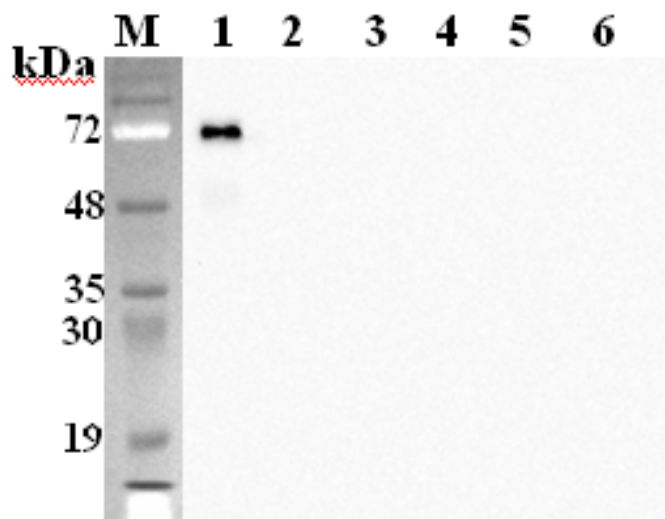


Figure: Western blot analysis of mouse Dlk1 using DLK1 (mouse), mAb (PF183E) (Prod. No. ALX-804-743) at 1:2'000 dilution.

1. DLK1 (mouse):Fc (human) (rec.) (Prod. No. ALX-201-416)
2. DLK1 (human):Fc (human) (rec.) (Prod. No. ALX-201-427)
3. DLL1 (mouse):Fc (human) (rec.) (Prod. No. ALX-201-455)
4. DLL4 (mouse):Fc (human) (rec.) (Prod. No. ALX-201-453)
5. Jagged-1 (mouse):Fc (human) (rec.) (Prod. No. ALX-201-463)
6. A unrelated protein (Fc fusion)

Handling & Storage

Use/Stability	Stable for up to 6 months when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Protein delta homolog 1, Pref-1, Preadipocyte factor 1, Fetal antigen 1, FA1
Application	ELISA, WB
Clone	PF183E
Crossreactivity	Does not cross-react with human DLK1.
Formulation	Liquid. 0.2µm-filtered solution in PBS, pH 7.4. Contains no preservatives.
Host	Rat
Immunogen	Recombinant mouse Dlk1 (extracellular domain).
Isotype	IgG2a
Recommendation Dilutions/Conditions	ELISA (direct and indirect, 1:2,000-1:10,000)Western Blot (1:2,000-1:10,000 using ECL; suggested blocking and dilution buffer is TBST containing 0.05% Tween 20 and 5% skim milk; suggested incubation time is 1 hour at room temperature)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse

UniProt ID

Q09163

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](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-
be@enzolifesciences.com](mailto:info-be@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)