DLK1 (mouse) monoclonal antibody (PF183E)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-743-C050

50µg

Manuals, SDS & CofA

View Online »

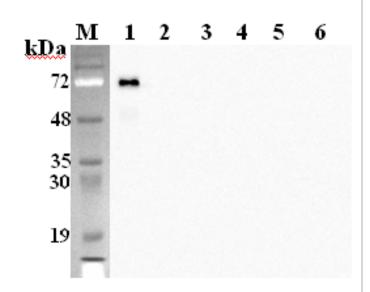


Figure: Western blot analysis of mouse Dlk1 using DLK1 (mouse), mAb (PF183E) (Prod. No. ALX-804-743) at 1:2'000 dilution.

- 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).

lgG2a

Recommendation Dilutions/ConditionsELISA (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

Q09163

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com