PPARγ2 (mouse) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

Ordering Information

Order Online »

ALX-210-192-R100

100µl

Manuals, SDS & CofA

View Online »

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 Peroxisome proliferator activated receptor y2

Application FUNC, Gel Supershift, ICC, IHC, WB

Application Notes Functional Application: Inhibits PPARy2 DNA binding

Detects a band of ~56kDa by Western blot.

Crossreactivity Does not cross-react with PPARα or PPARδ.

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 1-

> 16 (M1GETLGDSPIDPESDS16) of human PPARy2 (peroxisome proliferator activated receptor y2).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Immunocytochemistry (1:500)Immunohistochemistry

> (1:500)Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

UniProt ID P37231

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com