Guanylyl cyclase $\alpha_1^{}\beta_1^{}$ (soluble) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 12

View Online »

Ordering Information

Order Online »

ALX-210-786-R100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name sGCα1β1

Application WB

Application Notes Western Blot: Works on endogenous

Not recommended for the detection of human and mouse

 $sGC\alpha_1$ subunits, because of the lack of sensitivity.

Crossreactivity Does not cross-react with human and mouse $sGC\alpha 1$.

Formulation Liquid. In 10mM TRIS-HCl, pH 7.5, containing 150mM

sodium chloride and 0.02% thimerosal.

Host Chicken

Immunogen Purified rat lung soluble guanylyl cyclase α1β1

(sGCα1β1) (native protein).

Purity Detail Purified IgY from egg yolk.

Recommendation Dilutions/ConditionsWestern Blot(1:500)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

Specificity Recognizes endogenous rat $sGC\alpha_1\beta_1$. Recognizes human

and mouse $sGC\beta_1$.

UniProt ID P19686

Last modified: May 29, 2024



info-

eu@enzolifesciences.com