

Guanylyl cyclase $\alpha_1\beta_1$ (soluble) polyclonal antibody

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

Citations: 12

[View Online »](#)

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 sGC α 1 β 1

Application WB

Application Notes Western Blot: Works on endogenous
Not recommended for the detection of human and mouse sGC α ₁ subunits, because of the lack of sensitivity.

Crossreactivity Does not cross-react with human and mouse sGC α 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 α 1 β 1 (sGC α 1 β 1) (native protein).

Purity Detail Purified IgY from egg yolk.

Recommendation Dilutions/Conditions 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 Human, Mouse, Rat

Specificity Recognizes endogenous rat sGC α ₁ β ₁. Recognizes human and mouse sGC β ₁.

UniProt ID P19686

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