

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 sGCα ₁ subunits, because of the lack of sensitivity.
Crossreactivity	Does not cross-react with human and mouse sGCα1.
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/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 sGCα ₁ β ₁ . Recognizes human and mouse sGCβ ₁ .
UniProt ID	P19686



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 1 72 440 1335
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

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