

Uteroglobulin recombinant monoclonal antibody (122)

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

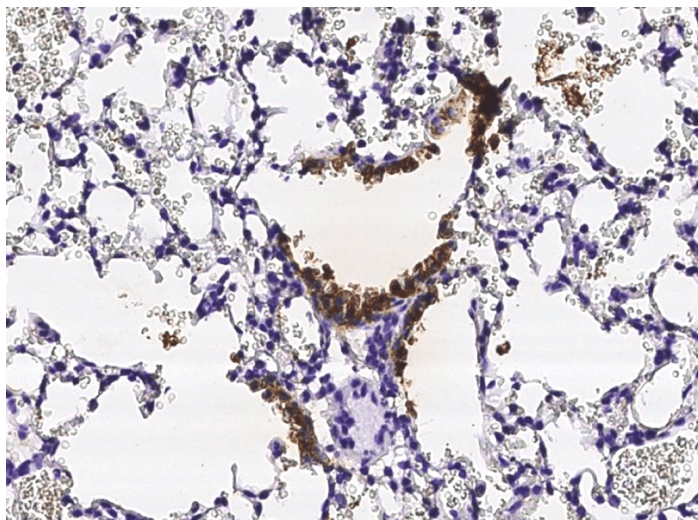
Ordering Information

[Order Online »](#)

ENZ-ABS1014-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of mouse SCGB1A1 in mouse lung with rabbit monoclonal antibody (1:200, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling	Protected from light. Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	SCGB1A1
Application	IHC (PS)
Clone	122
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant mouse Uteroglobin / SCGB1A1 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:100-1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: September 17, 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