

# Smooth muscle actin monoclonal antibody (CGA7) (Ready-to-Use)

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-30933	~6 µg - 6ml
-----------	-------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application IHC (PS), WB

Clone CGA7

Crossreactivity Does not cross-react with actin of skeletal or cardiac muscle cells.

Formulation Liquid. *Ready-to-use* solution in PBS containing 0.1% sodium azide and 1% BSA.

Host Mouse

Immunogen Partially purified chicken gizzard actin.

Isotype IgG2b

Purity Detail Affinity purified.

**Recommendation**  
**Dilutions/Conditions** Immunohistochemistry (paraffin sections; optimal results with this antibody are obtained when the tissue to be stained is fixed in alcohol-based fixatives such as methacarn. Aldehyde-fixed (formalin, e.g.) specimens are not recommended for use with CGA7. Choose a good quality enzyme-linked anti-mouse IgG2b antibody for detection) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Chicken, Human, Monkey, Rabbit

**Specificity** Recognizes actin isotypes  $\alpha$  and  $\gamma$  of smooth muscle cells. This antibody labels leiomyosarcomas and leiomyomas but does not react with other neoplasms such as carcinoma or lymphoma. Except for myoepithelium, it is nonreactive with other mesenchymal and all epithelial cells.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-bel@enzolifesciences.com](mailto:info-bel@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

**Worry-free Guarantee**

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