

Smooth muscle actin monoclonal antibody (CGA7)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-C34933	~10 µg - 0.5ml
------------	----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°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. 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).General suggested starting dilution is 1:20-1:50.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 α and γ 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.

Worry-free Guarantee

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



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
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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