

c-Fos polyclonal antibody

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-130-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, IHC (FS), IHC (PS), IP, WB

Formulation Liquid. In 0.05M sodium phosphate buffer containing 0.1% sodium azide and 0.2% gelatin.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 4-17(S4GFNADYEASSSRC17) of human c-Fos.

Purity Detail Immunoaffinity purified IgG.

Recommendation Dilutions/Conditions Immunocytochemistry (1-5µg/ml)Immunohistochemistry (frozen sections, 5-10µg/ml)Immunohistochemistry (paraffin sections, 5-10µg/ml, pepsin or heat pre-treatment required)Immunoprecipitation (1-2µg per reaction)Western Blot (1-5µg/ml).Suggested starting concentrations are provided.Antibodies should be titrated for optimal results in individual systems.

Species Reactivity Human, Mouse, Rat, Virus

Specificity Recognizes c-Fos and v-Fos (viral oncogene Fos).

UniProt ID P01100

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)