

c-Fos polyclonal antibody

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

Citations: 1

[View Online »](#)

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

Last modified: May 29, 2024

info@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420

info@enzolifesciences.com

France
Phone: +33 472 440 655

info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526

info@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488

Phone: +44 1392 825900

info@enzolifesciences.com