

CMV native antigen

Ordering Information

[Order Online »](#)

ALX-630-125-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Suitable for antibody capture
- High titre levels
- Offers low background interference

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Sonicate/homogenize briefly before use.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Cytomegalovirus native antigen
Appearance	Cloudy liquid.
Formulation	Liquid. In 0.2% Triton X-100, 10mM Tris-HCl, 1mM EDTA and 100mM NaCl, pH 7.8.
Purity	≥95%
Source	From cytomegalovirus, strain AD169. Inactivated by 0.2% Triton X-100.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)