

CLIP monoclonal antibody (CerCLIP.1)

CLIP (class II-associated invariant chain peptide) is an important peptide in the MHC-II antigen processing pathway, formed by the enzymatic cleavage of amino acids 81-104 from the invariant chain. During MHC-II synthesis in the ER and transport to endocytic vesicles, the antigen binding site of MHC-II is obstructed by the presence of the invariant chain. As the MHC-II complex moves through the Golgi, the invariant chain is degraded, leaving the CLIP fragment in the MHC-II peptide binding groove. CLIP is eventually displaced by the loading of antigenic peptides, followed by presentation on the APC cell surface to CD4+ T-cells.

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-CSA-600-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, IP, WB

Clone CerCLIP.1

Formulation Liquid. In PBS containing 0.09% sodium azide.

GenBank ID K01144

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of human CLIP.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Flow Cytometry (1:200) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from mouse ascites.

Species Reactivity Human

UniProt ID P04233

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)