

# 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 Dilutions/Conditions	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](#)

.



Last modified: May 29, 2024

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-be@enzolifesciences.com](mailto:info-be@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-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)