

# CD45 (human) clonal antibody (E19-G) (Ready-to-Use)

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

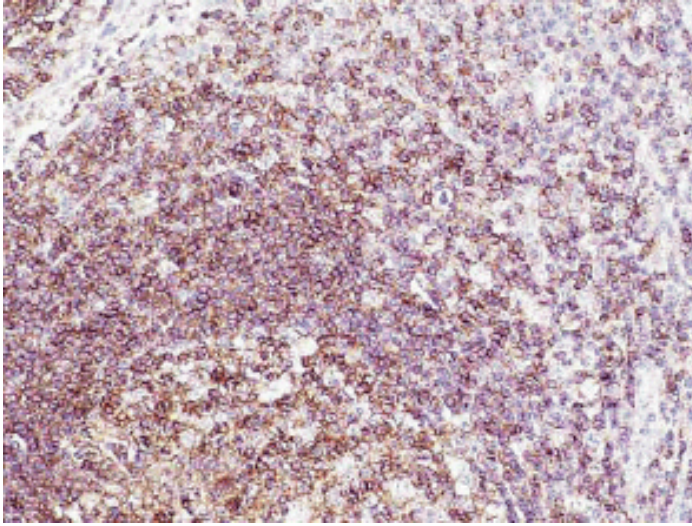
## Ordering Information

[Order Online »](#)

ALX-810-410-L015	15ml
------------------	------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Formalin-fixed and paraffin-embedded human tonsil tissue (4µm) stained with cAb to CD45 (human) (E19-G) shows strong positive immunostaining. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.

## Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored as recommended.
Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	IHC (PS)
Clone	E19-G
Formulation	Liquid. <i>Ready-to-use</i> solution in 20mM TRIS/HCl, pH 8.0, containing 20mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a C-terminal sequence of human CD45.
Isotype	IgG
Purity Detail	Major clone of IgGs obtained from immunoaffinity purified immunoglobulins corresponding to immunogenic peptide.
Species Reactivity	Human

## Technical Info / Product Notes

**Clonal Rabbit Antibody** is a pure homogenous fraction of rabbit immunoglobulin (IgG) which corresponds to a unique linear epitope on the target protein. This epitope is designed after a detailed analysis of the antigen structure, regulatory post-translational modifications, protein/protein interaction domain and protein folding. The chosen sequence (peptide) is synthesized and the rabbit is immunized. The clone is then separated from the crude antiserum by “*in vitro* cloning technology”. The resulting product is monospecific and characterized by superior specificity, affinity and avidity.

UniProt ID



Worry-free Guarantee

P08575

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  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-eu@enzolifesciences.com](mailto:info-eu@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)

This antibody is covered by our **Worry-Free Guarantee**.