

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 »](#)

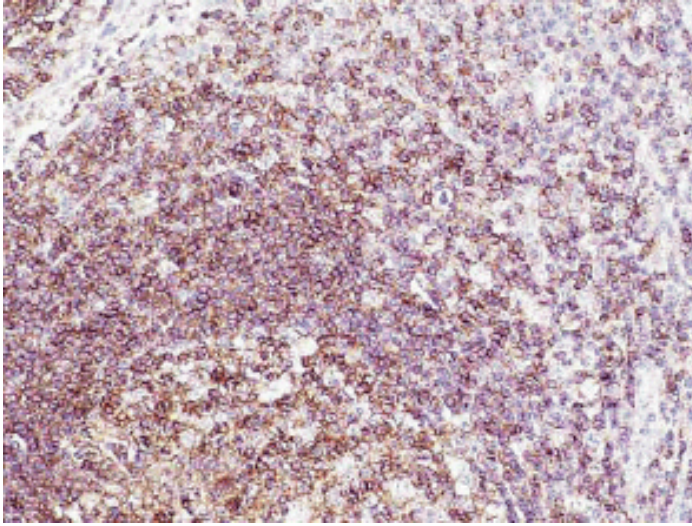


Figure: Formalin-fixed and paraffin-embedded human tonsil tissue (4µm) stained with cAb to CD45 (human) (E19-G) shows strong positive immunostaining. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.

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

P08575

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
[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)