

CD74 monoclonal antibody (LN2)

CD74, also known as the invariant chain, is involved in MHC class II antigen presentation. It is expressed on B cells and various epithelial tumors, including certain subtypes of breast and kidney cancers.

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

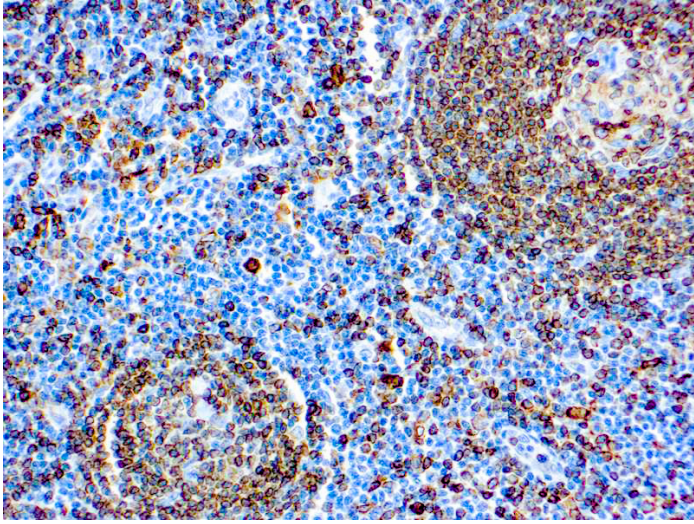
Ordering Information

[Order Online »](#)

ENZ-ABS1087-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human tonsil stained with CD74 antibody

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HLA class II histocompatibility antigen γ chain
Application	IHC (PS)
Application Notes	Pretreatment – EDTA Buffer pH 8.0. Incubation & Temperature 30 min @ RT
Clone	LN2
Formulation	Contains sodium azide as a preservative.
Host	Mouse
Immunogen	SU-DHL-4 lymphoma cells.
Isotype	IgG1k
Positive Control	Squamous Lung Carcinoma, Tonsil
Purity Detail	Purified immunoglobulin.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:50-1:200) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: June 6, 2025



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)