

CD6 recombinant monoclonal antibody (C6/2884R)

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

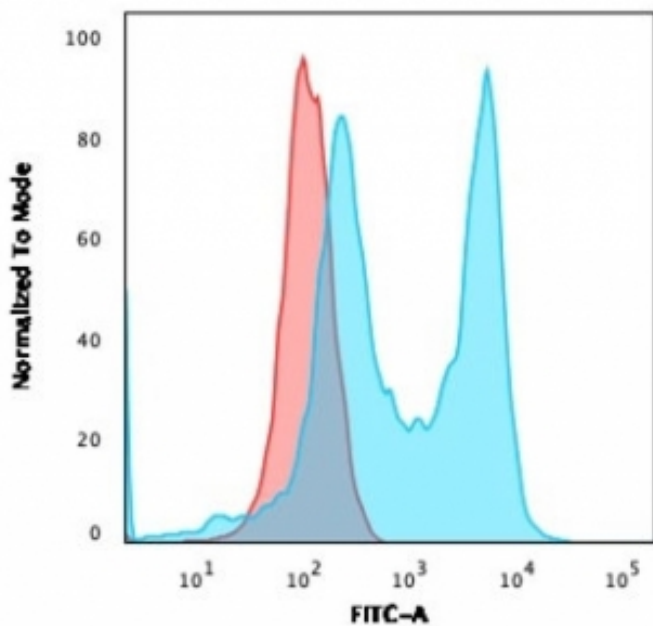
Ordering Information

[Order Online »](#)

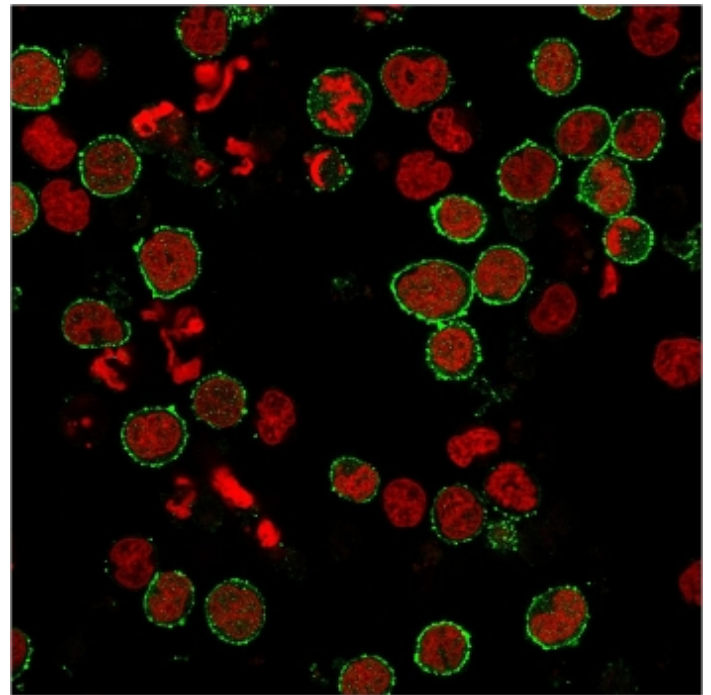
ENZ-ABS881-0100	100µg
-----------------	-------

Manuals, SDS & CofA

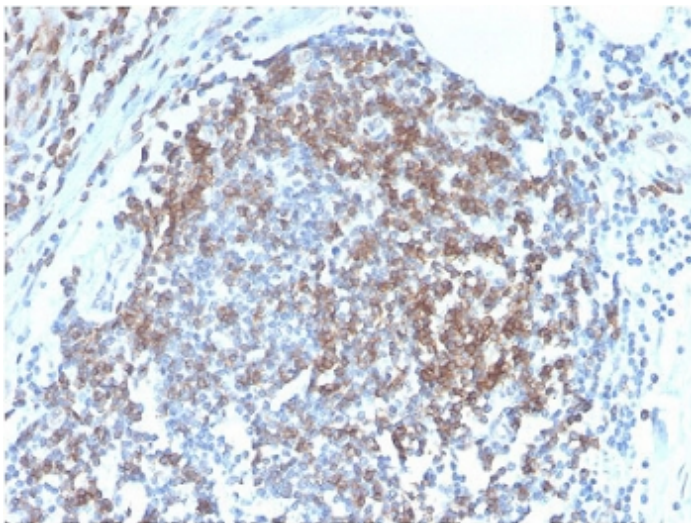
[View Online »](#)



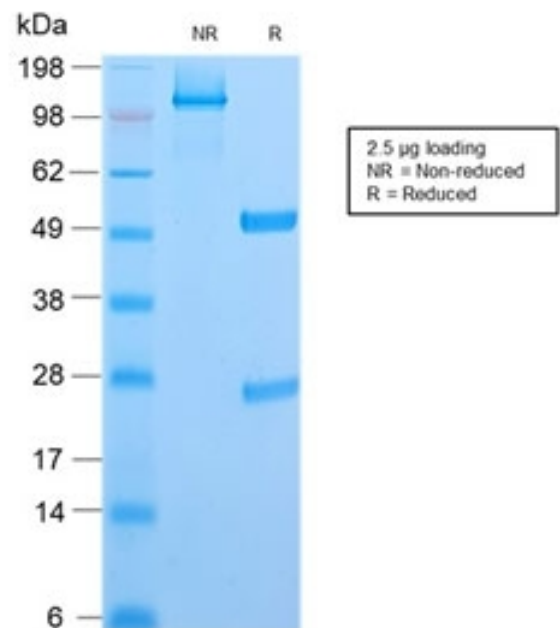
Flow cytometry testing of human MOLT4 cells with CD6 antibody (clone C6/2884R); Red=isotype control, Blue= CD6 antibody.



Immunofluorescent staining of human MOLT4 cells with CD6 antibody (clone C6/2884R, green) and Reddot nuclear stain (red).



IHC staining of FFPE human tonsil tissue with recombinant CD6 antibody (clone C6/2884R). HIER: boil tissue sections in pH6, 10mM citrate buffer, for 10-20 min followed by cooling at RT for 20 min.



SDS-PAGE analysis of purified, BSA-free recombinant CD6 antibody (clone C6/2884R) as confirmation of integrity and purity.

Handling & Storage

Handling	Store the recombinant CD6 antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide).
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Flow Cytometry, IF, IHC (PS)
Clone	C6/2884R
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Recombinant human CD6.
Isotype	IgGk
Purity Detail	Protein A affinity chromatography purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1-2µg/ml)Immunofluorescence (1-2µg/ml)Immunohistochemistry (FFPE) (0.5-1µg/ml for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P30203

Worry-free Guarantee

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

.



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

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

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

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

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

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

Last modified: July 12, 2024