

CD52 Recombinant monoclonal antibody (YTH 34.5-G2b) (Rat IgG2bκ)

Human CD52 antigen, also known as CAMPATH-1. CD52 is expressed at high density by lymphocytes, monocytes, eosinophils, thymocytes and macrophages. It is expressed by most lymphoid derived malignancies, although expression on myeloma cells is variable.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CDw52, Cambridge pathology 1 antigen, Epididymal secretory protein E5, Human epididymis-specific protein 5
Application	Cytotoxicity studies, ELISA, Flow Cytometry, IHC (FS), IHC (PS), in vivo Assay, WB
Clone	YTH 34.5-G2b (Campath-1G)
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rat
Immunogen	Human peripheral mononuclear cells enriched for T-cells.
Isotype	IgG2bk
Purity Detail	Protein A affinity purified.
Species Reactivity	Human, Rhesus monkey
UniProt ID	P31358

Worry-free Guarantee

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



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)

Last modified: May 29, 2024