

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](#).

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS455-0200 | 200µg |
|-----------------|-------|

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-
fr@enzolifesciences.com](mailto:info-fr@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