

CD19 Recombinant monoclonal antibody (HD37) (Mouse IgG1κ)

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

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS417-0200 | 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 | B-lymphocyte antigen CD19, B-lymphocyte surface antigen B4, Differentiation antigen CD19 T-cell surface antigen Leu-12 |
| Application | Flow Cytometry, FUNC, IF, IHC, IP |
| Clone | HD37 |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Mouse |
| Immunogen | Cells from a hairy cell leukemia patient. |
| Isotype | IgG1κ |
| Purity Detail | Protein A affinity purified. |
| Species Reactivity | Human |
| UniProt ID | P15391 |

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024