

# CD30 Recombinant monoclonal antibody (Ki-4) (Mouse IgG1κ)

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

|                             |  |
|-----------------------------|--|
| <b>Alternative Name</b>     | TNFRSF8, Tumor necrosis factor receptor superfamily member 8, CD30L receptor, Ki-1 antigen, Lymphocyte activation antigen CD30 |
| <b>Application</b>          | ELISA, Flow Cytometry, FUNC, IF, IHC   |
| <b>Application Notes</b>    | Functional Application: Inhibition   |
| <b>Clone</b>                | Ki-4   |
| <b>Formulation</b>          | Liquid. In PBS containing 0.02% Proclin 300.   |
| <b>Host</b>                 | Mouse  |
| <b>Immunogen</b>            | Human CD30.  |
| <b>Isotype</b>              | IgG1κ  |
| <b>Purity Detail</b>        | Protein A affinity purified.   |
| <b>Species Reactivity</b>   | Human  |
| <b>UniProt ID</b>           | P28908   |
| <b>Worry-free Guarantee</b> | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>   |

Last modified: May 29, 2024



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)