

# His(6) Recombinant monoclonal antibody (Rabbit IgGκ)

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

## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ENZ-ABS528-0200 | 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               | Polyhistidine tag                                                                                                                                                      |
| Application                    | ELISA, WB                                                                                                                                                              |
| Clone                          | 3D5                                                                                                                                                                    |
| Formulation                    | Liquid. In PBS containing 0.02% Proclin 300.                                                                                                                           |
| Host                           | Rabbit                                                                                                                                                                 |
| Immunogen                      | Purified his tagged Fab fragment.                                                                                                                                      |
| Isotype                        | IgGk                                                                                                                                                                   |
| Purity Detail                  | Protein A affinity purified.                                                                                                                                           |
| Specificity                    | Recognizes carboxy-terminal oligo-histidine tags (6x His tags) of fusion proteins. It also has affinity for internal His-tags but no affinity for N-terminal His-tags. |
| Technical Info / Product Notes | This is a chimeric rabbit antibody, based on the original mouse IgG2b format, created for improved compatibility with existing reagents, assays and techniques.        |
| Worry-free Guarantee           | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .                                                                                                 |



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)