

# CD14 (mouse), (recombinant) (His- tag)

The myeloid differentiation antigen CD14 acts as the major receptor for bacterial LPS. The dominant form of the recombinant wildtype CD14 ist the 50kDa protein.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-134-C010	10µg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	The reconstituted protein should be used immediately or stored in undiluted working aliquots at -20°C. For long term storage it is recommended to add a carrier protein (0.1% BSA).
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Biological Activity</b>	Inhibits binding of LPS to CD14. Up to 5µg/ml mouse CD14 inhibit binding of FITC-LPS (0.5µg/ml) to mouse CD14 <sup>+</sup> CHO transfectants (FACS).
<b>Concentration</b>	0.25mg/ml after reconstitution.
<b>Endotoxin Content</b>	<0.00015EU/µg protein (LAL test; Charles River).
<b>Formulation</b>	Lyophilized from PBS, pH 7.2.
<b>MW</b>	~45kDa
<b>Purity</b>	≥90-95% (SDS-PAGE)
<b>Purity Detail</b>	His-tag-affinity purified.
<b>Reconstitution</b>	Reconstitute with 40µl sterile, ultra-pure distilled water. This solution can then be diluted with PBS or other buffers.
<b>Source</b>	Produced in CHO cells.
<b>UniProt ID</b>	P10810



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)