

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 »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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).
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

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