CD14 (mouse), (recombinant) (Histag)

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

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

