

MRC1 (human), (recombinant) (His- tag)

Ordering Information

[Order Online »](#)

ENZ-PRT328-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa M

200
150
120
100
85
70
60
50
40
30
25
20



The recombinant human LOXL2 consists of 760 amino acids and has a predicted molecular mass of 85.5 kDa.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD206, Macrophage mannose receptor 1, MMR

Endotoxin Content <1 EU/μg (LAL method)

Formulation Lyophilized from sterile PBS, pH 7.4.

MW ~168.3 kDa (SDS-PAGE)

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in HEK293 cells. A DNA sequence encoding the human MRC1 (Leu19-Ser1386) was expressed with a polyhistidine tag at the C-terminus.

UniProt ID P22897

Last modified: November 7, 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)