

PRODUCT INFORMATION



CD163L1 SRCR8 Domain (human, recombinant)

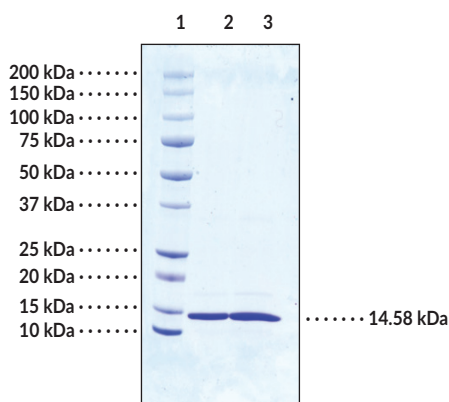
Item No. 41208

Overview and Properties

Synonyms:	CD163B, CD163L1, CD163 Antigen B, CD163 Antigen-like 1, CD163 Molecule-like 1, hCD163L SRCR8, M160, Scavenger Receptor Cysteine-rich Type 1 Protein M160
Source:	Recombinant human C-terminal His-tagged CD163L1 SRCR8 domain expressed in insect cells
Amino Acids:	790-895
Molecular Weight:	14.58 kDa
Storage:	-80°C (as supplied)
Stability:	≥6 months
Purity:	≥85% estimated by SDS-PAGE
Supplied in:	20 mM Tris, pH 8.0, 100 mM sodium chloride
Protein Concentration:	<i>batch specific</i> mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: CD163L1 SRCR8 Domain (2 µg)
Lane 3: CD163L1 SRCR8 Domain (4 µg)

SDS-PAGE Analysis of CD163L1 SRCR8 Domain.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/16/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

CD163 molecule-like 1 (CD163L1), also known as CD163 antigen B (CD163B), is an extracellular membrane receptor and a member of the scavenger receptor cysteine-rich (SRCR) family.^{1,2} It is composed of an N-terminal ectodomain, which contains 12 disulfide-rich SRCR class B domains that mediate protein-protein interactions, a transmembrane segment, and a short cytoplasmic tail.¹ Alternative splicing of CD163L1 mRNA produces two variants that differ in length of the cytoplasmic tail.^{1,3} CD163L1 is found in monocytes and M2 macrophages and can be induced by colony-stimulating factor 1 (CSF1) and IL-10 in macrophages.³ Replacement of porcine CD163L1 SRCR5 domain with human CD163L1 SRCR8 domain increases resistance to porcine reproductive and respiratory syndrome virus 1 (PRRSV1) infection in pigs.⁴ Cayman's CD163L1 SRCR8 Domain (human, recombinant) protein consists of 105 amino acids and has a calculated molecular weight of 14.58 kDa.

References

1. Grønlund, J., Vitved, L., Lausen, M., *et al.* Cloning of a novel scavenger receptor cysteine-rich type I transmembrane molecule (M160) expressed by human macrophages. *J. Immunol.* **165**(11), 6406-6415 (2000).
2. Etzerodt, A. and Moestrup, S.K. CD163 and inflammation: Biological, diagnostic, and therapeutic aspects. *Antioxid. Redox Signal.* **18**(17), 2352-2363 (2013).
3. González-Domínguez, É., Samaniego, R., Flores-Sevilla, J.L., *et al.* CD163L1 and CLEC5A discriminate subsets of human resident and inflammatory macrophages in vivo. *J. Leukoc. Biol.* **98**(4), 453-466 (2015).
4. Wells, K.D., Bardot, R., Whitworth, K.M., *et al.* Replacement of porcine CD163 scavenger receptor cysteine-rich domain 5 with a CD163-like homolog confers resistance of pigs to genotype 1 but not genotype 2 porcine reproductive and respiratory syndrome virus. *J. Virol.* **91**(2), e01521-e015316 (2017).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM