

PRODUCT INFORMATION



CD20 Extracellular Domain (human, recombinant; aa 141-188)

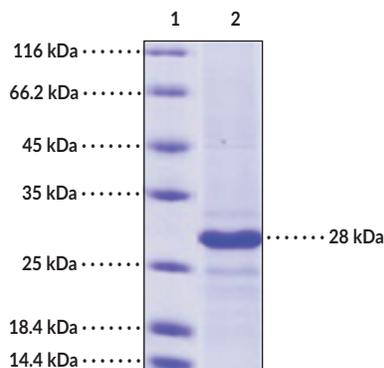
Item No. 44202

Overview and Properties

Synonyms:	B Lymphocyte Cell Surface Antigen B1, Membrane-spanning 4-domains Subfamily A Member 1
Source:	Recombinant human N-terminal TrxA-tagged CD20 extracellular domain (141-188) expressed in <i>E. coli</i>
Amino Acids:	141-188
Uniprot No.:	P11836
Molecular Weight:	23.9 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥80% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 50 mM Tris, 1mM EDTA, and 5% glycerol, pH 8.0
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: CD20 Extracellular Domain

SDS-PAGE Analysis of CD20 Extracellular Domain. This protein has a calculated molecular weight of 23.9 kDa. It has an apparent molecular weight of approximately 28 kDa by SDS-PAGE under reducing conditions.

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, 01/06/2026

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

CD20 is a non-glycosylated protein encoded by *MS4A1* in humans.¹ It is composed of four transmembrane domains, a single intracellular loop, and two extracellular loop domains with both the N and C-termini located in the cytosol. CD20 is a general B cell marker that is expressed from the late pre-B lymphocyte stage but is not expressed by pro-B lymphocytes and is lost in terminally differentiated plasma cells and plasmablasts. It forms supramolecular complexes with CD53, CD81, and CD82, as well as MHCII, CD40, B cell receptors, and C-terminal Src kinase-binding protein (CBP) to contribute to signal transduction. *MS4A1* expression is variable in B cell malignancies, with the lowest expression found in patients with chronic lymphocytic leukemia (CLL) and the highest expression found in patients with diffuse large B cell lymphoma (DLBCL) or hairy cell lymphomas. *MS4A1* expression is enriched on IFN- γ -inducible T-box transcription factor-expressing B cells in blood isolated from patients with multiple sclerosis.² Cayman's CD20 Extracellular Domain (human, recombinant; aa 141-188) protein consists of 218 amino acids and has a calculated molecular weight of 23.9 kDa. By SDS page, under reducing conditions, the apparent molecular mass of the protein is approximately 28 kDa.

References

1. Pavlasova, G., and Mraz, M. The regulation and function of CD20: An "enigma" of B-cell biology and targeted therapy. *Haematologica* **105(6)**, 1494-1506 (2020).
2. van Langelaar, J., Rijvers, L., Smolders, J., *et al.* B and T cells driving multiple sclerosis: Identity, mechanisms and potential triggers. *Front. Immunol.* **11**, 760 (2020).

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