

PRODUCT INFORMATION



2B4/CD244 Extracellular Domain (human, recombinant)

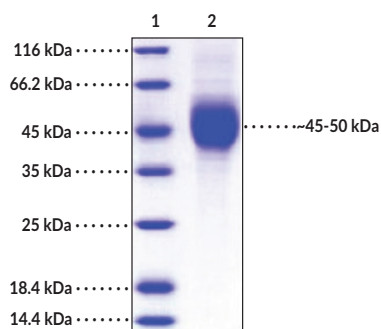
Item No. 31820

Overview and Properties

Synonyms:	Cluster of Differentiation 244, NAIL, Natural Killer Cell Receptor 2B4, NK Cell Activation-inducing Ligand, NKR2B4, Nmrk, SLAM Family Member 4, SLAMF4
Source:	Active recombinant human C-terminal His-tagged 2B4 expressed in HEK293 cells
Amino Acids:	22-221
Uniprot No.:	Q9BZW8
Molecular Weight:	23.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥95% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Bioactivity:	See figure for details

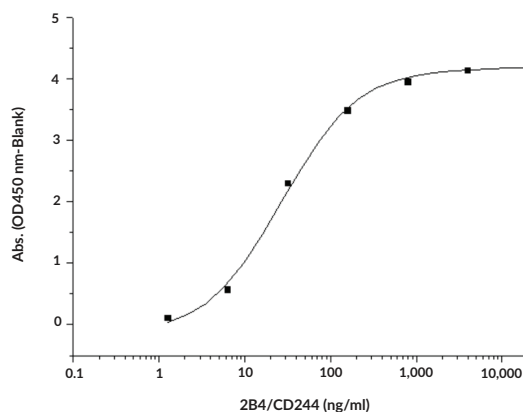
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: 2B4/CD244

SDS-PAGE Analysis of 2B4/CD244. This protein has a calculated molecular weight of 23.8 kDa. It has an apparent molecular weight of approximately 45 to 50 kDa by SDS-PAGE under reducing conditions due to apparent glycosylation.



2B4/CD244 Functional ELISA. Immobilized human 2B4/CD244 Extracellular Domain (human, recombinant) at 2 μg/ml (100 μl/well) can bind human CD48. The EC₅₀ value of human CD48 is 0.39 μg/ml.

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, 10/16/2020

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

PRODUCT INFORMATION



Description

2B4, also known as CD244, is a cell surface receptor and a member of the signaling lymphocytic activation molecule (SLAM) receptor family (SLAMF).¹ It is comprised of an extracellular domain containing immunoglobulin (Ig) variable-like (IgV) and Ig constant 2-like (IgC2) motifs and several glycosylation sites, a membrane-spanning region, and a cytoplasmic domain containing serine, tyrosine, and threonine phosphorylation sites.² It is expressed on natural killer (NK) cells, certain T cells, monocytes, basophils, and eosinophils. 2B4 has an activating role when the adapter protein SLAM-associated protein (SAP) is present and bound to intracellular tyrosine-based switch motifs (ITSM) on the cytoplasmic domain.¹ When SAP is absent, SHP-1, SHP-2, SHIP-1, or CsK are able to bind, which leads to inhibitory signaling. 2B4 has *cis*- and *trans*-interactions with the cell surface receptor CD48 during the initiation and continuation of allergic reactions. It has a complex role in cancer, with male, but not female, *CD244*^{-/-} mice rejecting tumor cells in a xenograft model, however, its role is primarily inhibitory in tumor-associated immune cells.² The expression of *CD244* is decreased in NK cells from patients with multiple myeloma or systemic lupus erythematosus (SLE).¹ 2B4 is associated with X-linked lymphoproliferative disease (XLPD), which is characterized by a mutation in SAP that prevents it from binding to 2B4.³ Cayman's 2B4/CD244 Extracellular Domain (human, recombinant) protein can be used for ELISA. This protein consists of 211 amino acids, has a calculated molecular weight of 23.8 kDa, and a predicted N-terminus of Cys22 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is approximately 45-50 kDa due to glycosylation.

References

1. Pahima, H., Puzzovio, P.G., and Levi-Schaffer, F. 2B4 and CD48: A powerful couple of the immune system. *Clin. Immunol.* **204**, 64-68 (2019).
2. Buller, C.W., Mathew, P.A., and Mathew, S.O. Roles of NK cell receptors 2B4 (CD244), CS1 (CD319), and LLT1 (CLEC2D) in cancer. *Cancers (Basel)* **12**(7), 1755 (2020).
3. Mathew, P.A. Regulation of NK cell function by 2B4 (CD244) receptor. *Trends Immunol.* **9**, 25-32 (2008).

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