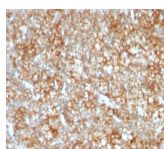


CD147 Monoclonal Mouse Antibody (CB43)



Product Description

CD147 is a transmembrane glycoprotein of the immunoglobulin superfamily. It is expressed more intensely on thymocytes than on mature peripheral blood T cells. CD147 is important in spermatogenesis, embryo implantation, neural network formation, and tumor progression. CD147 is involved in the regulation of matrix remodeling at the epidermal-dermal interface. It stimulates the production of interstitial collagenase, gelatinase A, stromelysin-1 and various metalloproteinases (MMPs) by fibroblasts. These enzymes, which are typically increased during tissue degradation and wound healing, are important factors in cancer invasion and metastasis.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Catalog number key for antibody number 0363, Anti-CD147 (CB43)

Product attributes

Antibody number	#0363
Antibody reactivity (target)	CD147
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CB43
Isotype	IgM
Molecular weight	35 kDa (non-reduced); 40 kDa (reduced)
Synonyms	5A11 Antigen; 5F7; Basigin; Blood brain barrier HT7 antigen BSG; Extracellular Matrix Metalloproteinase Inducer (EMMPRIN); Leukocyte Activation Antigen M6; Neurothelin; OK Blood Group Antigen; TCSF; Tumor Cell-derived Collagenase Stimulatory Factor
Human gene symbol	BSG
Entrez gene ID	682
SwissProt	P35613
Unigene	501293
Immunogen	Native human leucocytes
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	HSB2 cells. Renal Cell, Ovarian carcinoma or Melanoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Cancer, Developmental biology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	CF@405S Features
BNC88	CF@488A	490/515	488	GFP, FITC	CF@488A Features
BNC68	CF@568	562/583	532, 561	RFP, TRITC	CF@568 Features
BNC94	CF@594	593/614	561	Texas Red®	CF@594 Features
BNC40	CF@640R	642/662	633-640	Cy@5	CF@640R Features
BNC47	CF@647	650/665	633-640	Cy@5	CF@647 Features
BNC74	CF@740	742/767	633-685	775/50	CF@740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 29, 2025 at 01:40:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd147-cb43/>