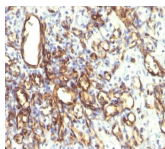


Podocalyxin Monoclonal Mouse Antibody (2A4)



Product Description

Podocalyxin is a member of the CD34 transmembrane sialomucin family. It is highly on kidney podocyte foot projections and plays essential roles in kidney development and homeostasis, blood filtration and urine formation. It is also expressed on vascular endothelia, hematopoietic progenitors and a subset of neurons. Overexpression of podocalyxin may be linked to more aggressive tumor behavior. Podocalyxin antibody can identify podocytes in the urine (podocyturia) that may indicate glomerular disease, pre-eclampsia, and other kidney pathology.

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 0483, Anti-Podocalyxin (2A4)

Product attributes

Antibody number	#0483
Antibody reactivity (target)	Podocalyxin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	2A4
Isotype	IgM
Molecular weight	165-170 kDa
Synonyms	PODXL, PDXL, GCTM-2; Gp200; PCLP1; Pcx; Podocalyxin-like protein 1
Human gene symbol	PODXL
Entrez gene ID	5420
SwissProt	O00592
Unigene	732423
Immunogen	A recombinant protein fragment containing the intracellular, transmembrane, and part of the extracellular domain of human podocalyxin
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human, Rabbit, Rat
Antibody application notes	Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	HeLa, Raji, Jurkat cells, Angiosarcoma, Breast, Prostate, Liver, Pancreatic, Kidney
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	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.
Cell/tissue expression	Kidney

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 December 16, 2025 at 10:10:44 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-podocalyxin-2a4/>