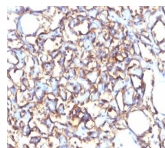


Podocalyxin Monoclonal Mouse Antibody (3D3)



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. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0427, Anti-Podocalyxin (3D3)**

Product attributes

Antibody number	#0427
Antibody reactivity (target)	Podocalyxin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3D3
Isotype	IgG1
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
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)
Species reactivity	Human, Rabbit, Rat
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Immunohistochemistry 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. Western blotting 0.5-1 ug/mL. Immunoprecipitation 1-2 ug/500 ug protein. 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. Protect fluorescent conjugates from light. 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
Cell/tissue expression	Kidney

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

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, LI-COR Bioscience.