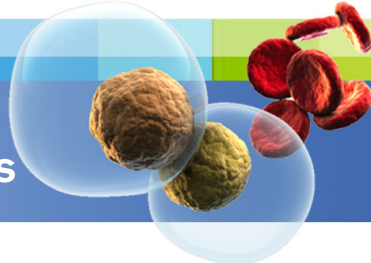




SINGLE-COLOR ANTIBODIES



NEW PRODUCT

CD138-Krome Orange

PN B42019

Antibody

Specificity: Human
Clone: B-A38
Isotype: IgG1 mouse

Reagent

Line: IOTest
Format: Liquid
Size: 0.5 mL
Status: ASR

Antigen characteristics

Other name: Syndecan-1,
Heparan sulfate proteoglycan

Gene Location: 2p24.1
Gene Name: SDC1
Molecular Weight: 30.5 kDa

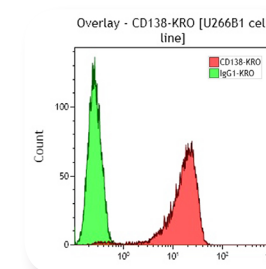
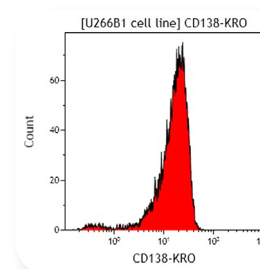
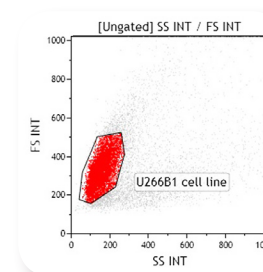
CD138 / syndecan-1 protein backbone is a single chain molecule. Five putative GAG attachment sites exist in the extracellular domain. The addition of GAG chains to the protein backbone of syndecans is constitutive and critical, since these provide all of the known extracellular ligand binding sites on syndecans. In general, GAG fine structure appears to reflect the cellular source of the syndecan. Expression of CD138 / syndecan-1 in human hematopoietic cells is restricted to plasma cells in normal bone marrow. Peripheral blood monocytes, T- and B-lymphocytes are not reactive to CD138. Tonsil cells and early B-cell precursors in human bone marrow also are CD138-negative. CD138 / syndecan-1 is also expressed in a variety of mature tissues, including epithelial cells, endothelial cells, fibroblasts, keratinocytes and normal hepatocytes. CD138 / syndecan-1 is proposed to have a variety of roles depending on the expressing cell types, and on the size and composition of carried heparan sulfates. It may act as an essential regulator of ligand-specific activation of primary signalling receptors at the cell surface, by co-engagement of various accessory ligands.

Formats available

	Size	Line	Status	Part #		Size	Line	Status	Part #
PE	100 tests	IOTest	CE	A54190	APC	0.5 mL	IOTest	ASR	A87787
PC5	100 tests	IOTest	CE	A54191	Pacific Blue*	0.5 mL	IOTest	ASR	B37788
PC5.5	0.5 mL	IOTest	ASR	A98835	Krome Orange	0.5 mL	IOTest	ASR	B42019

ASR: Analyte Specific Reagent. Analytical and performance characteristics are not established.
CE: CE-marked product

Example of results



U266B1 transfected cell line

References

- Carey, D., «Syndecans: Multifunctional cell-surface co-receptors», 1997, Biochem. J., 327, 1-16.
- Wijdenes, J., Clement, C., Klein, B., Dore, J.M., «CD138 (syndecan-1) Workshop panel report», 1997, Leucocyte Typing VI, White Cell Differentiation Antigens. Kishimoto, T., et al, Eds., Garland Publishing, Inc., 249-252.
- Subramanian, S.V., Fitzgerald, M.L., Bernfield, M., «Regulated shedding of syndecan-1 and -4 ectodomains by thrombin and growth factor receptor activation», 1997, J. Biol. Chem., 272, 14713-14720.



Life Sciences