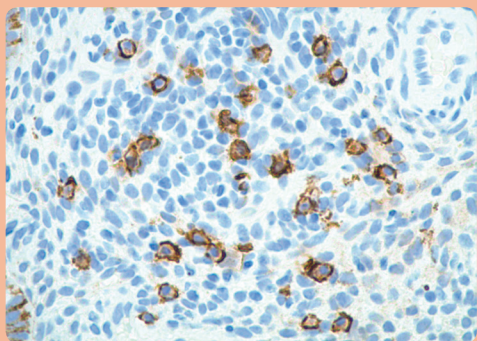


CD138 Monoclonal Antibody (ENZ-ABS683)

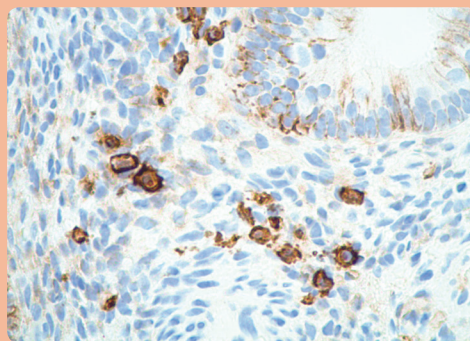
Cluster of differentiation 138 (CD138), also known as Syndecan-1, is a transmembrane glycoprotein present on the surface of cells during late stage differentiation. Anti-CD138 is used to differentiate marginal zone lymphoma from lymphoplasmacytic lymphoma. ALK+ Large B-Cell Lymphoma (LBCL) cells commonly stain positively for CD138, but not for CD20 and CD79a. Anti-CD138 reacts positively with HHV8-associated primary effusion lymphoma-expressing cells that lack B-cell markers. CD138 is also a useful marker for identifying and enumerating benign, reactive, or malignant plasma cells from bone marrow biopsies.

Enzo's CD138 mAb (ENZ-ABS683)



Endometrial Tissue – IHC staining was performed on an automated Leica Bond III

Competitor's CD138 mAb (mouse)



Endometrial Tissue – IHC staining was performed on an automated Leica Bond III

RELATED PRODUCTS

Product Name	Product #	Product Name	Product #
POLYVIEW® PLUS AP (Anti-Rabbit) Reagent	ENZ-ACC110	HIGHDEF® Blue IHC Chromogen (HRP)	ADI-950-151
POLYVIEW® PLUS AP (Anti-Mouse) Reagent	ENZ-ACC114	HIGHDEF® Green AP Chromogen/Substrate	ENZ-ACC130
POLYVIEW® PLUS HRP (Anti-Mouse) Reagent	ENZ-ACC104	HIGHDEF® Hematoxylin	ENZ-ACC106
POLYVIEW® PLUS HRP (Anti-Rabbit) Reagent	ENZ-ACC103	HIGHDEF® Red IHC Chromogen (AP)	ADI-950-140/1
SAVIEW® PLUS AP Reagent	ENZ-ACC111	HIGHDEF® Yellow IHC Chromogen (HRP)	ADI-950-170
SAVIEW® PLUS HRP Reagent	ENZ-ACC102	IHC/ISH Peroxidase Block	ENZ-ACC107
HIGHDEF® DAB Chromogen/Substrate Set	ENZ-ACC105	Antigen Retrieval Reagent, pH 6 (10X)	ENZ-ACC112
HIGHDEF® AEC (Red) IHC Chromogen (HRP)	ADI-950-210	Antigen Retrieval Reagent, pH 9 (10X)	ENZ-ACC113
HIGHDEF® Black IHC Chromogen (HRP)	ADI-950-171	Antibody Blocker/Diluent	ENZ-ACC108
HIGHDEF® Blue IHC Chromogen (AP)	ADI-950-150	Streptavidin Blocker/Diluent	ENZ-ACC109

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com

For Research Use Only
Not for Use in Diagnostic Procedures

© 2018 Enzo Life Sciences.