



# PRODUCT DATA SHEET

## Caldesmon Recombinant monoclonal antibody (RMCDN1-1)

ENZ-ABS336

### Product Number/Sizes

ENZ-ABS336-0100 100 µg

### Product Details

ALTERNATIVE NAME:	CAD, CDM
CLONE:	RMCDN1-1
HOST:	Rabbit
ISOTYPE:	IgGκ
IMMUNOGEN:	Recombinant full length human Caldesmon.
UNIPROT ID:	Q05682
SPECIES REACTIVITY:	Human
APPLICATIONS:	IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS:	Immunofluorescence (0.5-1 µg/mL) Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL for 30 min at RT) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
PURITY DETAIL:	Protein A affinity purified.
FORMULATION:	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING:	Blue Ice
LONG TERM STORAGE:	+4°C
SCIENTIFIC BACKGROUND:	Caldesmon is an actin- and myosin-binding protein implicated in the regulation of actomyosin interactions in smooth muscle and nonmuscle cells (could act as a bridge between myosin and actin filaments). Stimulates actin binding of tropomyosin which increases the stabilization of actin filament structure. In muscle tissues, inhibits the actomyosin ATPase by binding to F-actin. This inhibition is attenuated by calcium-calmodulin and is potentiated by tropomyosin. Interacts with actin, myosin, two molecules of tropomyosin and with calmodulin. Also play an essential role during cellular mitosis and receptor capping. Involved in Schwann cell migration during peripheral nerve regeneration.
REGULATORY STATUS:	RUO - Research Use Only

Revised 19-Dec-23

#### GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.  
10 Executive Blvd  
Farmingdale, NY 11735  
USA  
T 1-800-942-0430  
T 1-631-694-7070  
F 1-610-941-9252  
E [info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)

#### EUROPE/ASIA

Enzo Life Sciences (ELS) AG  
Industriestrasse 17, Postfach  
CH-4415 Lausen  
Switzerland  
T +41/061 926 89 89  
T +41/061 926 89 79  
E [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)