



PRODUCT DATA SHEET

Calponin Recombinant monoclonal antibody (RMCN1-1)

ENZ-ABS341

Product Number/Sizes

ENZ-ABS341-0100 100 µg

Product Details

CLONE: RMCN1-1
HOST: Rabbit
ISOTYPE: IgGκ
IMMUNOGEN: Recombinant human calponin.
UNIPROT ID: P51911
SPECIES REACTIVITY: Human
APPLICATIONS: Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10⁶ cells)
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: Calponin is a calmodulin, F-actin and tropomyosin binding protein which is thought to be involved in the regulation of smooth muscle contraction.
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
www.enzolifesciences.com

EUROPE/ASIA

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