



# PRODUCT DATA SHEET

## Calponin Recombinant monoclonal antibody (CNN1-1408R)

ENZ-ABS350

### Product Number/Sizes

ENZ-ABS350-0100 100 µg

### Product Details

CLONE: CNN1/1408R  
HOST: Rabbit  
ISOTYPE: IgGκ  
IMMUNOGEN: Recombinant full length human calponin-1.  
UNIPROT ID: P51911  
SPECIES REACTIVITY: Human, Rat  
APPLICATIONS: Flow Cytometry, IF, IHC (PS)  
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10<sup>6</sup> cells)  
Immunofluorescence (0.5-1 µg/mL)  
Immunohistochemistry (formalin-fixed paraffin sections) (1-2 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 20-Jul-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  
F +41/061 926 89 79  
E [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)