
CAPN2 Polyclonal Antibody, HRP Conjugated

(Catalog # A52891)

Background

Calcium-regulated non-lysosomal thiol-protease which catalyze limited proteolysis of substrates involved in cytoskeletal remodeling and signal transduction.

Description

CAPN2 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P17655

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Calpain-2 catalytic subunit protein (20-700AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Calpain M-type, Calpain large polypeptide L2, Calpain-2 large subunit, Millimolar-calpain, CAPN2, CANPL2

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
