
ARMC8 Polyclonal Antibody

(Catalog # A58154)

Background

Component of the CTLH E3 ubiquitin-protein ligase complex that selectively accepts ubiquitin from UBE2H and mediates ubiquitination and subsequent proteasomal degradation of the transcription factor HBP1.

Description

ARMC8 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q8IUR7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Armadillo repeat-containing protein 8 protein (1-240AA)

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

Armadillo repeat-containing protein 8, ARMC8, S863-2

Application

ELISA
