
MED27 Polyclonal Antibody

(Catalog #A65690)

Background

Component of the Mediator complex, a coactivator involved in the regulated transcription of nearly all RNA polymerase II-dependent genes. Mediator functions as a bridge to convey information from gene-specific regulatory proteins to the basal RNA polymerase II transcription machinery. Mediator is recruited to promoters by direct interactions with regulatory proteins and serves as a scaffold for the assembly of a functional preinitiation complex with RNA polymerase II and the general transcription factors.

Description

MED27 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

Q6P2C8

Purification

Affinity purified

Immunogen

Recombinant Human Mediator of RNA polymerase II transcription subunit 27 protein (39-108AA)

Storage

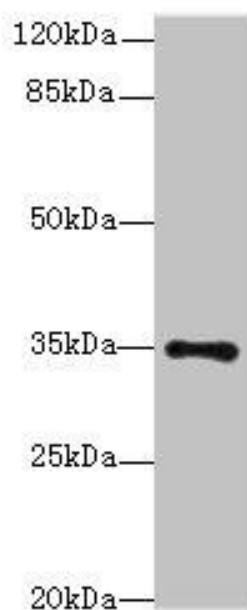
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Mediator of RNA polymerase II transcription subunit 27, Cofactor required for Sp1 transcriptional activation subunit 8, CRSP complex subunit 8, Mediator complex subunit 27, P37 TRAP/SMCC/PC2 subunit, Transcriptional coactivator CRSP34, MED27, CRSP34

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: MED27 Polyclonal Antibody at 0.5µg/ml + A431

whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 36, 32, 15 kDa

Observed band size: 36 kDa