

HSPA6 Polyclonal Antibody

(Catalog #A53804)

Description

HSPA6 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

P17066

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Heat shock 70 kDa protein 6 protein (1-637AA)

Storage

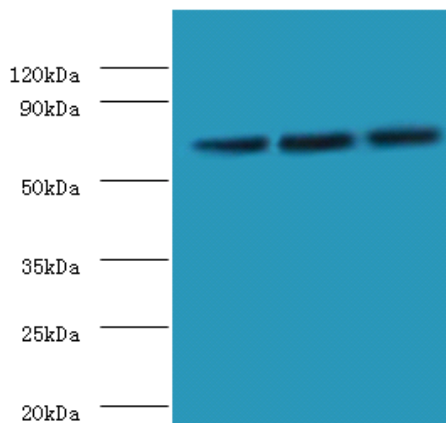
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Heat shock 70 kDa protein B' HSPA6 HSP70B'

Application

ELISA, WB; Recommended dilution: ELISA 1:500-1:1000, WB 1:1000 - 1:5000



Western blot

All lanes: HSPA6 Polyclonal Antibody at 2ug/ml

Lane 1: MCF-7 whole cell lysate

Lane 2: A431 whole cell lysate

Lane 3: Jurkat whole cell lysate

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 71kDa

Observed band size: 71kDa