

ANTI-PROTEASOME 20S SUBUNIT BETA 10

RABBIT MONOCLONAL ANTIBODY



CATALOG #: 3322

RECOMBINANT ANTIBODY

ALIASES

PSMB10; Proteasome 20S Subunit Beta 10; LMP10; MECL1; Proteasome (Prosome, Macropain) Subunit, Beta Type, 10; Multicatalytic Endopeptidase Complex Subunit MECL-1; Proteasome Subunit Beta Type-10; Low Molecular Mass Protein 10; Proteasome Subunit Beta 10; Proteasome Subunit Beta-2i; Macropain Subunit MECL-1; Proteasome MECL-1; EC 3.4.25.1; MGC1665; Beta2i; Proteasome Catalytic Subunit 2i; Proteasome Subunit Beta 7i; Proteasome Subunit Beta2i; Proteasome Subunit MECL1; Proteasome Subunit B2i; PRAAS5; BETA2I

BACKGROUND

Gene Name: *PSMB10*

UniProt Entry: P40306

NCBI Gene Entry: 5699

APPLICATION INFORMATION

Molecular Weight: Predicted, 29 kDa, observed, 25 kDa

Clonality: Rabbit monoclonal antibody

Clone ID: 24GB5050

Species Reactivity: Human

Applications Tested: Western blotting (WB)

IMMUNOGEN

A synthesized peptide derived from human PSMB10

SUPPORT

SUPPORT@GENUINBIOTECH.COM

TEL: +1-540-855-7041

ORDERS

SALES@GENUINBIOTECH.COM

FAX: +1-540-855-7041

WWW.GENUINBIOTECH.COM

ANTI-PROTEASOME 20S SUBUNIT BETA 10

RABBIT MONOCLONAL ANTIBODY

PAGE 2

ISOTYPE

Rabbit IgG

STORAGE BUFFER

Supplied in PBS (pH 7.4) containing 50% glycerol, and 0.02% sodium azide.

STORAGE

Store at -20 °C for one year.

RECOMMENDED DILUTIONS

Western Blotting (WB): 1:1,000–1:5,000

PROTOCOLS

For general and specific antibody protocols please visit our website. Read all instructions before using this product.

Note: This product is for research use only.

SUPPORT

SUPPORT@GENUINBIOTECH.COM

TEL: +1-540-855-7041

ORDERS

SALES@GENUINBIOTECH.COM

FAX: +1-540-855-7041

WWW.GENUINBIOTECH.COM

ANTIBODY VALIDATION

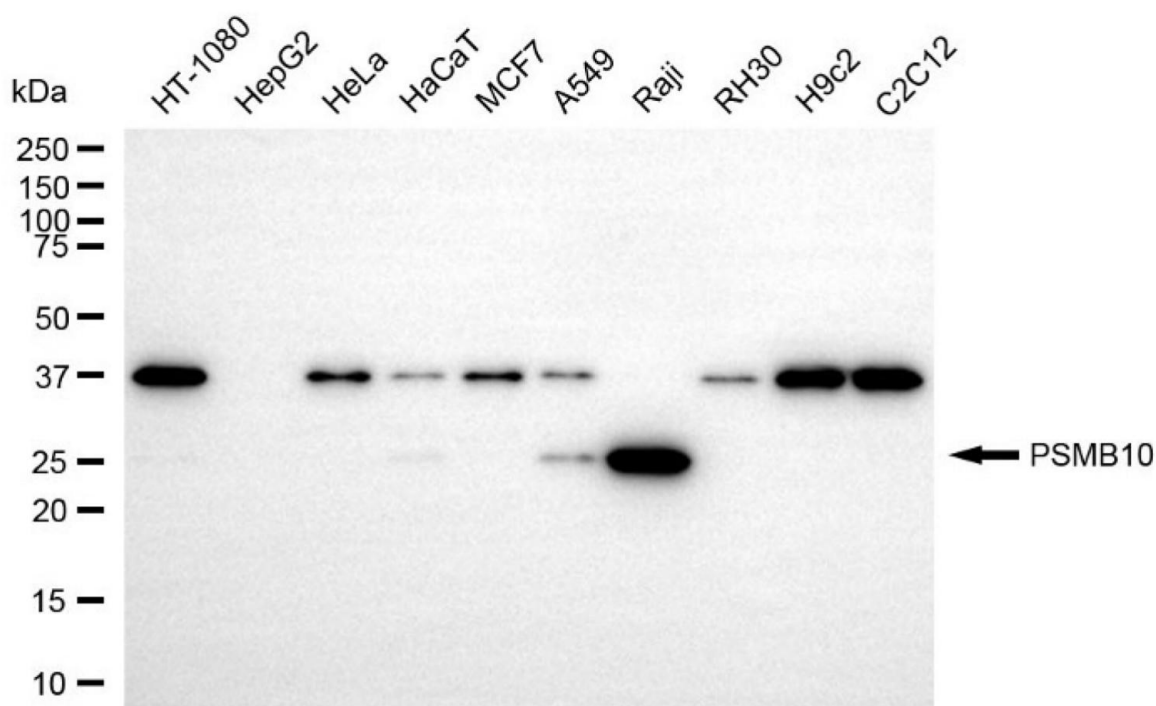


Figure 1. Western blotting analysis using anti-PSMB10 antibody (Cat#3322). Total cell lysates (30 μ g) from various cell lines were loaded and separated by SDS-PAGE. The blot was incubated with anti-PSMB10 antibody (Cat#3322, 1:5,000) and HRP-conjugated goat anti-rabbit secondary antibody (Cat#201, 1:20,000) respectively. Image was developed using NaQ™ ECL Substrate Kit (Cat#716). PSMB10, proteasome 20S subunit beta 10.

SUPPORT

SUPPORT@GENUINBIOTECH.COM

TEL: +1-540-855-7041

ORDERS

SALES@GENUINBIOTECH.COM

FAX: +1-540-855-7041

WWW.GENUINBIOTECH.COM