POMP polyclonal antibody

The assembly of individual mammalian proteasome subunits into catalytically active 20S proteasome is a complex process, with biogenesis of the mammalian 20S proteasome occuring \emph{via} precursor complexes containing α and unprocessed β subunits. The pro-peptides of the β subunits are thought to prevent premature conversion of the precursor complexes into matured particles and are needed for efficient β subunit incorporation. The complex biogenesis is tightly regulated, requiring additional components such as the maturation factor POMP (proteasome maturation protein), an ubiquitous protein in eukaryotic cells. Mammalian POMP is functionally related to but not interchangable with its yeast homologue Ump1. POMP is associated purely with the precursor intermediates and is degraded upon final maturation. Northern blot experiments have revealed that expression of POMP is induced after treatment with interferon γ .

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-PW9715-0025	25μΙ
BML-PW9715-0100	100μΙ

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome maturation protein, UMP1 homolog

Application IHC, IP, WB

Formulation Liquid. In PBS containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 129-141 of human

POMP.

Purity Detail Purified by salt precipitation and peptide

immunochromatography.

Species Reactivity Human, Monkey, Rat

UniProt ID Q9Y244

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024

