

---

# Proteasome 20S Alpha6 Subunit Monoclonal Antibody [MCP20]

(Catalog No. A-8425)

---

**Description:**

Mouse monoclonal antibody to Proteasome 20S alpha6 subunit, Clone MCP20

**Size:**

100 µl

**Formulation**

Liquid. In PBS containing 10mM sodium azide.

**Immunogen**

Human placenta derived proteasomes

**Specificity:**

Human, Rabbit

**Isotype:**

IgG1

**Purification:**

Partially purified ascites

**Storage:**

Store at -20°C. Avoid multiple freeze/thaw cycles.

**Application:**

IHC, IP, WB

**Ordering Information****Products**

Proteasome 20S Alpha6 Subunit Monoclonal Antibody [MCP20]

**Size**

100 µl

**Cat. No.**

A-8425-100

*This product is for research purposes only. Not intended for use in diagnostic procedures.*