

## HP1 $\gamma$ antibody (mAb)

---

**Catalog No:** 39981

**Clone:** 2MOD-1G6

**Isotype:** IgG1k

**Application(s):** ICC, IF, WB

**Reactivity:** Human, Mouse

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 0.57  $\mu\text{g}/\mu\text{l}$

**Quantity:** 100  $\mu\text{g}$

**Molecular Weight:** 21 kDa

---

**Background:** Heterochromatin Protein 1 gamma (HP1 gamma) is one of the mammalian homologues of Drosophila HP1, first identified through its role in the maintenance of regions of the genome in a heterochromatic state. HP1 gamma contains a chromo domain and a chromo shadow domain. Through the chromo domain, HP1 proteins are recruited to regions of chromatin that exhibit methylation of lysine 9 of histone H3.

**Immunogen:** This HP1 $\gamma$  antibody was raised against recombinant full-length mouse HP1 $\gamma$ .

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

Published Applications:

ICC/IF

WB

See reference for experimental details. Individual optimization may be required.

**References:** HP1 $\gamma$  antibody has been cited in one or more publications. To see an up-to-date listing of papers that describe its use, please go to [www.activemotif.com/catalog/details/39981](http://www.activemotif.com/catalog/details/39981).

**Storage and Guarantee:** Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.