

## CAF-1 p150 antibody (mAb)

---

**Catalog Nos:** 39997, 39998

**RRID:** AB\_2793424

**Clone:** 1

**Isotype:** IgG1

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 150 kDa

---

**Background:** CAF-1 p150 (Chromatin assembly factor-1 p150 subunit, CAF1A, CHAF1A) is the largest subunit of the CAF-1 complex that is required for the assembly of nucleosomes onto newly replicated DNA. Specifically, CAF-1 assists in the loading of the H3-H4 tetramer and may be involved in DNA damage repair.

**Immunogen:** The CAF-1 p150 antibody was raised against full-length CAF-1 purified from human 293 cells.

**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:

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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