

PCNA antibody (mAb)

Catalog Nos: 61079, 61080

RRID: AB_2793499

Clone: 16D10

Isotype: IgG2b

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

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 32 kDa

Background: PCNA (Proliferating cell nuclear antigen) is a marker for cells in early G1 phase and S phase of the cell cycle. It is found in the nucleus and is a cofactor of DNA polymerase delta. PCNA acts as a homotrimer and helps increase the processivity of leading strand synthesis during DNA replication.

Immunogen: This PCNA antibody was raised against full-length recombinant human PCNA.

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:

IP

ICC/IF

WB

ELISA

See references for more information. Individual optimization may be required.

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.