

PRODUCT INFORMATION



PCNA Monoclonal Antibody (Clone IPO-38)

Item No. 10004805

Overview and Properties

Contents:	This vial contains 100 µg of affinity-purified monoclonal antibody.
Synonyms:	Nuclear Antigen of Proliferating Cells, Proliferating Cell Nuclear Antigen
Species Reactivity:	(+) Multiple species PCNA (species- and tissue-unspecific)
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	200 µl PBS, pH 7.2, with 0.02% sodium azide
Clone:	IPO-38
Host:	Mouse
Isotype:	IgM
Applications:	Flow cytometry (FC), Immunocytochemistry (ICC), Immunohistochemistry (IHC) (frozen and paraffin-embedded sections), and Western blot (WB); the recommended starting concentration for WB is 1-2 µg/ml. FC, ICC, IHC, and other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

The proliferating cell nuclear antigen (PCNA) becomes detectable in the G₁ phase of the cell cycle and reaches a maximum level of expression during mitosis. PCNA is highly conserved across species and is a critical part of the DNA polymerase δ holoenzyme.¹⁻³ The level of PCNA expression depends on the proliferative potential of the cell examined. Cayman's PCNA Monoclonal Antibody can be used for analysis of PCNA by Western blot, flow cytometry, immunocytochemistry, and immunohistochemistry (frozen and paraffin-embedded tissue) on samples from multiple species. This monoclonal antibody stains proliferating cells in a wide range of normal tissues, as well as in tissues undergoing pathological processes and after mitogenic and allogeneic stimulations.

References

1. Tan, C.-K., Castillo, C., So, A.G., *et al.* An auxiliary protein for DNA polymerase-δ from fetal calf thymus. *J. Biol. Chem.* **261(26)**, 12310-12316 (1986).
2. Fukuda, K., Morioka, H., Imajou, S., *et al.* Structure-function relationship of the eukaryotic DNA replication factor, proliferating cell nuclear antigen. *J. Biol. Chem.* **270(38)**, 22527-22534 (1995).
3. Travali, S., Ku, D.-H., Rizzo, M.G., *et al.* Structure of the human gene for the proliferating cell nuclear antigen. *J. Biol. Chem.* **264(13)**, 7466-7472 (1989).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/11/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM