

PCAF Protein

Recombinant protein expressed in E. coli cells

Catalog # P09-31G

Lot # U014-1

Product Description

Recombinant human PCAF (431-end) was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [NM_003884](#).

Gene Aliases

CAF, GCN5, GCN5L, P/CAF, GCN5L1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

PCAF is a protein that shows homology to the yeast GCN5 and associates with p300/CBP. PCAF has in vitro and in vivo binding activity with CBP and p300, and competes with E1A for binding sites in p300/CBP (1). PCAF encodes a predicted 832-amino acid protein that showed by RNA blotting to be expressed in all tissues, most strongly in heart and skeletal muscle. PCAF has histone acetyl transferase activity with core histones and nucleosome core particles, indicating that this protein plays a direct role in transcriptional regulation. PCAF is required for MyoD activity and muscle differentiation. PCAF directly acetylated MyoD which then displays an increased affinity for its DNA target (2).

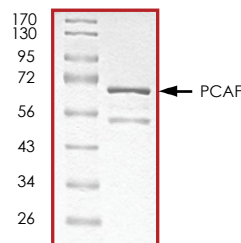
References

1. Yang, X J. et al: A p300/CBP-associated factor that competes with the adenoviral oncoprotein E1A. Nature 382: 319-324, 1996.
2. Sartorelli, V. et al: Acetylation of MyoD directed by PCAF is necessary for the execution of the muscle program. Molec. Cell 4: 725-734, 1999.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

Purity



The purity of PCAF was determined to be **>70%** by densitometry. Approx. MW **69kDa**.

PCAF Protein

Recombinant protein expressed in E. coli cells

Catalog Number **P09-31G**

Specific Lot Number **U014-1**

Purity **>70%**

Concentration **0.2 µg/µl**

Stability **1 yr at -70°C from date of shipment**

Storage & Shipping **Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.**