

---

## Pcaf Protein

(Catalog No. E29000)

---

### Description

Recombinant P/CAF encoding human sequence from 492 to 658. Genebank access no. Q92831)

### Molecular Weight (kDa)

19.3 kDa

### Formulation

Liquid. In 20mM sodium citrate, pH 6.0, containing 0.25M NaCl, 1mg/ml BSA and 10% glycerol.

### Source

E. Coli

### Size

50 µg

### Storage Temp & Conditions

-80°C; Centrifuge vial prior to opening

### Purification Notes

Purity : ≥95%

### Application Notes

Use in bioassay, or store at -80°C for future use.

### Ordering Information:

Products	Size	Cat. No.
Pcaf Protein	50 µg	E29000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*