

---

# PCBP1 Polyclonal Antibody

(Catalog # A72648)

---

## Background

This intronless gene is thought to have been generated by retrotransposition of a fully processed PCBP-2 mRNA. This gene and PCBP-2 have paralogues (PCBP3 and PCBP4) which are thought to have arisen as a result of duplication events of entire genes. The protein encoded by this gene appears to be multifunctional. It along with PCBP-2 and hnRNP-K corresponds to the major cellular poly(rC)-binding protein. It contains three K-homologous (KH) domains which may be involved in RNA binding. This encoded protein together with PCBP-2 also functions as translational coactivators of poliovirus RNA via a sequence-specific interaction with stem-loop IV of the IRES and promote poliovirus RNA replication by binding to its 5'-terminal cloverleaf structure. It has also been implicated in translational control of the 15-lipoxygenase mRNA, human Papillomavirus type 16 L2 mRNA, and hepatitis A virus RNA. The encoded protein is also suggested to play a part in formation of a sequence-specific alpha-globin mRNP complex which is associated with alpha-globin mRNA stability.

## Description

PCBP1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

## Specificity

Human, Mouse, Rat

## Isotype

IgG

## Uniprot ID

Q15365

## Purification

Affinity Purification

## Immunogen

A synthetic peptide corresponding to a sequence within amino acids 100-200 of human PCBP1 (NP\_006187.2).

## Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

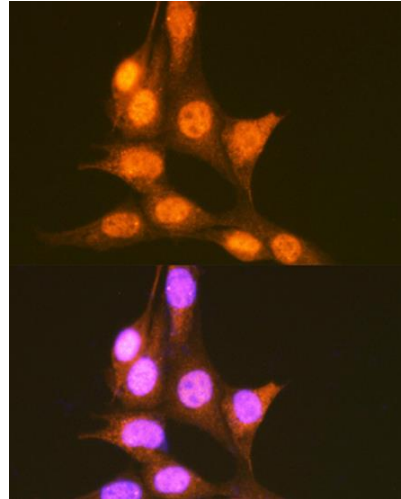
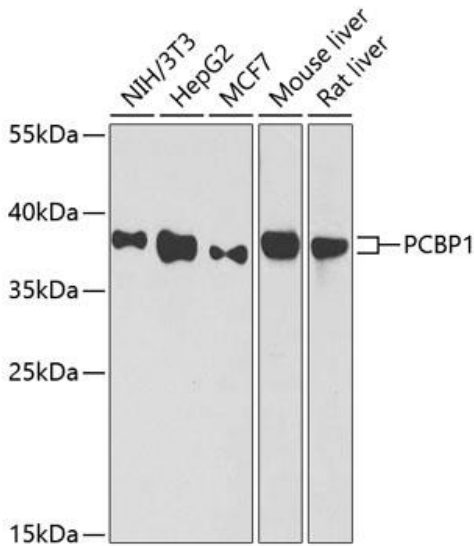
## Alternative Names

PCBP1; HEL-S-85; HNRPE1; HNRPX; hnRNP-E1; hnRNP-X; poly(rC)-binding protein 1

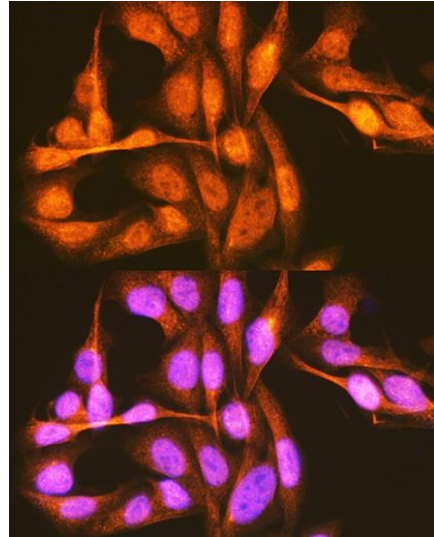
## Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100, IF 1:50 - 1:200

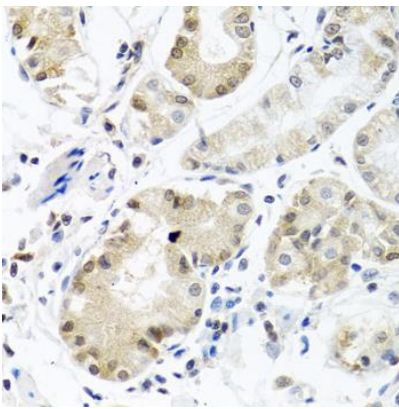
---



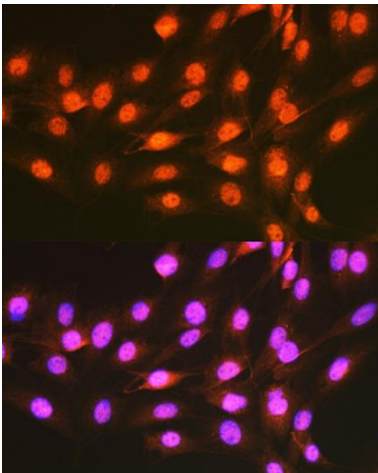
Immunofluorescence analysis of NIH-3T3 cells using PCBP1 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunofluorescence analysis of U-2 OS cells using PCBP1 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded human stomach using PCBP1 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunofluorescence analysis of C6 cells using PCBP1 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.