

## Recombinant CREB Protein

**Catalog No:** 31107

**Source:** Human

**Expressed In:** *E. coli*

**Contents:** 25 µg of Recombinant CREB Protein

### Protein Details

Recombinant human active CREB (amino acids 1-327) is produced in *E. coli* as a maltose binding protein (MBP) fusion protein with an amino terminal polyhistidine tag and purified via sequential chromatography to >95%. Recombinant CREB is phosphorylatable in vitro using either recombinant Protein Kinase A or Rsk immunoprecipitated from stimulated cells.

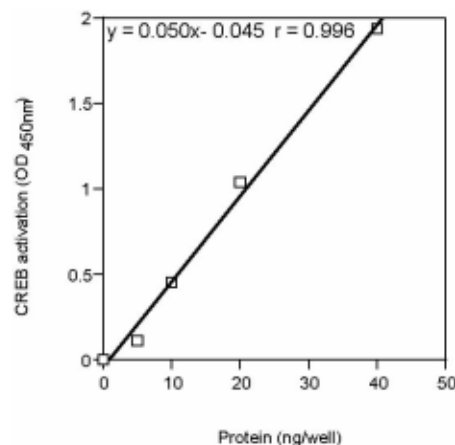
### Dilution of Protein

Briefly centrifuge the vial prior to opening to bring contents to the bottom of the tube. Reconstitute in PBS to a concentration of 0.2-1.0 mg/ml. Following reconstitution, store in aliquots at -80°C. If you are performing TransAM assays, dilute the recombinant protein using the Complete Lysis Buffer provided in the TransAM Kit.

### Application

Recombinant CREB is suitable for use in Western blotting and TransAM assays. The standard curve for TransAM™ CREB was generated using the range of 40-10 ng of protein. The molecular weight of CREB is ~81 kDa.

**NOTE:** The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.



The standard curve for TransAM was generated using a range of 40-10 ng of protein and run on the TransAM CREB ELISA Kit.

### Storage

Lyophilized proteins can be stored at 4°C, preferably desiccated. Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.

### Research use

This product is for research use only and is not for use in diagnostic procedures.