

Data Sheet

CREBBP (1081-1197), His-Tag

Human, recombinant, N-terminal His- tag

Catalog #: 31119

Lot#: 180622-1 **Conc.:** 1.36 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 10%
glycerol.

Stability: At least 6 months at -80°C

References:

1. Panagopoulos, I., *et al.*, *Hum Mol Genet.* 2001 Feb 15; **10(4)**: 395-404.
2. Huang, W.C., *et al.*, *Mol Cell.* 2007 Apr 13; **26(1)**:75-87.

Description:

Human bromodomain CREB binding protein (CREBBP), also known as CBP, RSTS and KAT3A, GenBank Accession No. NM_004380, a.a. 1081-1197, with N-terminal His-tag, MW=15 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE

Coomassie staining

Lane 1:

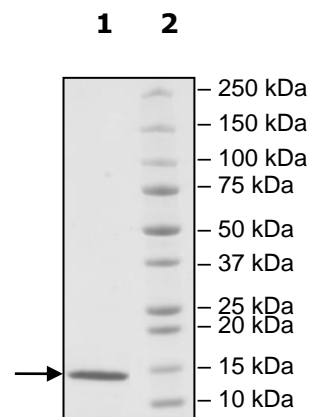
3 μg

Lane 2:

Protein Marker

MW: 15 kDa

Purity: 90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com