

CREB binding protein (catalytic domain) (human), (recombinant)

Although originally termed histone acetyltransferases (HATs) for their lysine acetylation activity on histone N-terminal tails, CBP and its paralogue, p300, have been shown to acetylate a variety of non-histone proteins including p53, DNA polymerase β and nuclear import factors. p300/CBP acetylations play regulatory roles in transcription, DNA repair and replication, the cell cycle, p53 turnover, and nuclear import. In addition to functions which overlap with those of p300, CBP has a number of unique functions, including a role in p27Kip1 induction.

Citations: 5

[View Online »](#)

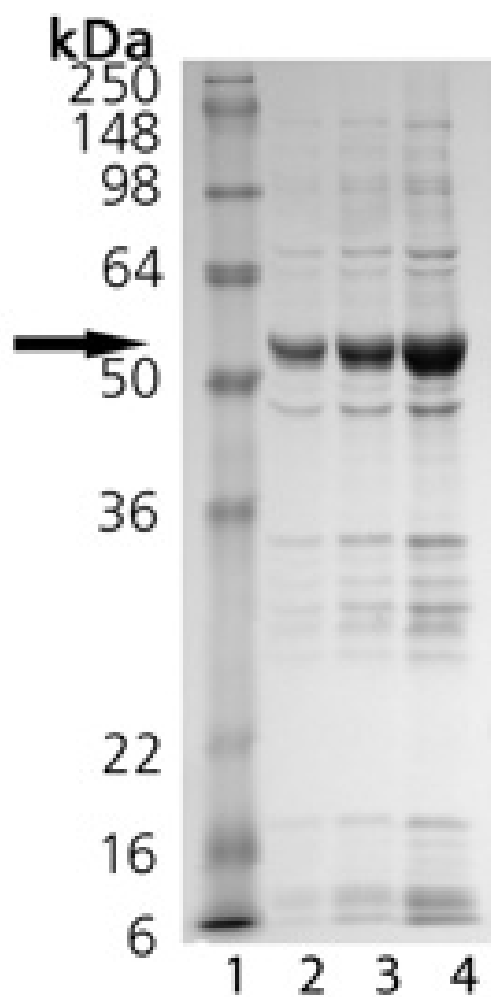
Ordering Information

[Order Online »](#)

BML-SE452-0100	100 μ g
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 µg, Lane 3: 2 µg, Lane 4: 5 µg CBP.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CBP
Application Notes	Can be used to study CBP kinetics and regulation, for screening for inhibitors or activators and for radiolabeling of proteins or peptides with, for example [³ H]-Acetyl CoA.
Formulation	Liquid. In 50mM TRIS-HCl, pH 8.0, containing 0.1mM EDTA, 10% glycerol.
MW	45.4 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in <i>E. coli</i> . Catalytic domain (aa 1319-1710) of human CBP.
Specific Activity	≥250 pmol/min/μg assayed as production of CoA-SH from AcCoA in the presence of a peptide comprising human p53 residues 368-386 (Prod. No. BML-P198). CoA-SH is determined colorimetrically by reaction with DTNB (5,5'-Dithiobis(2-nitrobenzoic acid)).
UniProt ID	Q92793

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)