

## Data Sheet

### ACC1, FLAG-His-tags

Human, recombinant, C-terminal FLAG-His-tags

**Catalog #:** 50202

**Lot #:** 170208

**Conc.:** 0.18 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0. 110 mM NaCl, 2.2 mM KCl, 0.1 mg/mL FLAG peptide and 20% Glycerol

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Sinilnikova, O.M. *et al.*, *Cancer Epidemiol. Biomarkers Prev.* **16 (3)**, 409-415 (2007)
2. Olsen, J.V. *et al.*, *Cell* **127 (3)**, 635-648 (2006)

**Description:** Human ACC1 (acetyl-Coenzyme A carboxylase alpha transcript variant 1) (GenBank Accession No. NM\_198834), amino acids 1–2383(end) with C-terminal FLAG and His-tags, MW = 271 kDa, expressed in Baculovirus infected Sf9 cell expression system.

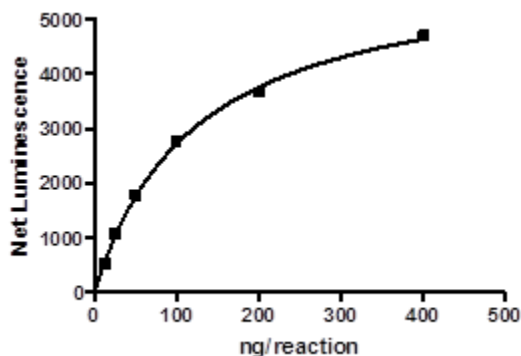
**Specific Activity:**  $\geq 58$  pmol/min/ $\mu\text{g}$ .

Assay was done using ADP-Glo (Promega) reagents with 20  $\mu\text{M}$  ATP, 20  $\mu\text{M}$  acetyl-CoA and 30 mM  $\text{NaHCO}_3$  in 25  $\mu\text{L}$  ACC assay buffer (30 mM HEPES pH=7.4, 2 mM  $\text{MgCl}_2$ , 1 mM DTT, and 2 mM Sodium citrate). Reaction was performed at room temperature for 40 min.

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

## Quality Assurance

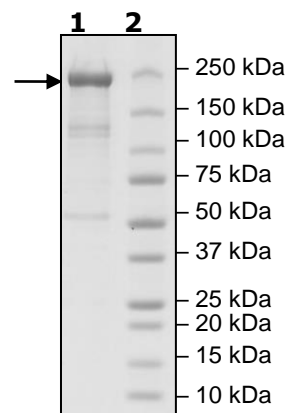
**Specific Activity**



**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
2  $\mu\text{g}$  ACC1  
**Lane 2:**  
Protein Marker

**MW:** 271 kDa  
**Purity:**  $\geq 67\%$



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court West, Suite B  
 San Diego, CA 92121  
 Tel: 1.858.829.3082  
 Fax: 1.858.481.8694  
 Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Amino Acid Sequence:**

MWWSTLMSILRARSFWKWISTQTVRIIRAVRAHFGGIMDEPSPLAQPLELNQHSRFIIGSVSEDNS  
 EDEISNLVKLLDLEEKESLSPASVGSDTLSDLGISLQDGLALHIRSSMSGHLHLVKQGRDRKKID  
 SQRDFTVASPAEFVTRFGGNKVIEKVLIANNGIAAVKCMRSIRRWSYEMFRNERAIRFVVMVTPED  
 LKANA EYIKMADHYVVPVPGPNNNNYANVELILDIAKRIPVQAVWAGWGHASENPKLPELLLNKGI  
 AFMGPPSQAMWALGDKIASSIVAQTAGIPTLPWSGSGLRVDWQENDFSKRILNVPQELYEKGYVKD  
 VDDGLQAAEEVGYPMVIKASEGGGKIRKVNADDPNLFRQVQAEVPGSPIFVMRLAKQSRHLE  
 VQILADQYGN AISLFG RDCSVQRRHQKIEEAPATIATPAVFEHMEQCAVKLAKMVGYSAGTVEY  
 LYSQDGSFYFLELNPRLQVEHPCTEMVADVNLPAAQLOIAMGIPLYRIKDIRMMYGVSPWGDSPID  
 FEDSAHVPCPRGHVIAARITSENPDGFKPSSGTVQELNFRSNKNVWGYFSVAAAGGLHEFADSQF  
 GHCFSWGENREEAISNMVVALKEL SIRGDFRTTVEYLIKLETESFQMNRI DTGWLDRLIAEKVQA  
 ERPDTMLGVVCGALHVADVSLRNSVSNFLHSLERGOVLP AHTLLNTVDVELIYEGVKYVLKVTRQS  
 PNSYVVMNGSCVEVDVHRLSDGGLLSYDGSSYTYMKEEVD RYRITIGNKTCVFEKENDP SVMR  
 SPSAGKLIQYIVEDGGHVFAGQCYAEIEVMKMVMTLTAVESGCIHYVKRPGAALDPGCVLAKMQLD  
 NPSKVQQAELHTGSLPRIQSTALRGEKLRHFVHYVLDNLVNMNGYCLPDPFFSSKVKDWVERLMK  
 TLRDPSLPLELQDIMTSVSGRIPPNVEKSIKKEMAQYASNITSVLCQFP SQQIANILDSHAATLN  
 RKSEREVFFMNTQSIVQLVQRYSGIRGHMKAVVMDLLRQYLRVETQFQNGHYDKCVFALREENKS  
 DMNTV LNYIFSHAQVTKNLLVTMLIDQLCGRDPTLTDELLN ILTELTQLSKTTNAKVALRARQVL  
 IASHLPSYELRHNVESIFLSAIDMYGHQFCIENLQKLI LSETSIFDVLNPFYHSNQVVRMAALE  
 VYVRRAYIAYELNSVQHRQLKDNTCVVEFQFMLPTSHPNRGNIP TLRMSFSSNLNHYGMTHVASV  
 SDVLLDNSFTPPCQRMGMVSRFTFEDFVRI FDEVMGCFSDSPPQSPTFPEAGHTSLYDEDK VPRD  
 EPIHILNVAIKTDCDIEDDRLAAMFREFTQQNKATLV DHGIRRLTFLVAQKDFRKQVNYEVDRRFH  
 REFPKFFTFRRARDKFEEDRIYRHLEPALAFQLELNRMRNFDLTAIPC ANHKMHLYLGA AKVEVGTE  
 VTDYRFFVRAIIRHSDLVTKEASFEYLQNEGERL LLEAMDELEVA FNNTNVRTDCNHI FLNFVPTV  
 IMDPSKIEESVRSVMRYGSRWLKLRVLQAE LKINIRLTPTGKAIPIRLFLT NESGYLDISLYKE  
 VTDSRTAQIMFQAYGDKQGPLHGMLINTPYVT KDLLQSKRFQAQSLGTTYIYDIPEMFRQSLIKLW  
 ESMSTQAF LPSPLPSDMLTYTELVLDDQGQLVHMNRLPGGNEIGMVAWKMTFKSPEYPEG RDIIV  
 IGNDITYRIGSFGPQEDLLFLRASELARAEGIPRIYVSANSGARIGLAE EIRHMFHVAWVDPEDPY  
 KGYRYLYLTPQDYKRVSALNSVHCEHVEDEGESRYKITDIIGKEEGIGPENLRGSGMIAGESSL AY  
 NEIITISLVT CRAIGIGAYLVRLGQRTIQVENS HLLITGAGALNKVLGREVYTSNNQLGGIQIMHN  
 NGVTHCTVCDDFEGVFTVLHWSYMPKSVHSSVPL LNSKDPIDRIIEFVPTKTPYDPRWMLAGRPH  
 PTQKGQWLSGFFDYGSFSEIMQPWAQTVVVGRARLGGIPVGVVAVETRTVELSIPADPANLDSEAK  
 IIQQAGQVWFPSAFKTYQAIKDFNREGLPLMV FANWRGFGSGMKDMDQVLKFGAYIVDGLRECC  
 QPVLVYI PPQAE LRGGSWVIDSSINPRHMEMYADRESRGSVLEPEGTVEIKFRKDLVKTMRRVD  
 PVYIHLAERLGTPELSTAERKELENK LKEREEFLIPIYHQVAVQFADLHDT PGRMQEKGVISDILD  
 WKTSRTFFYWRRLRRLLEDLVKKKIHNANPELTDGQIQAMLRRWFVEVEGTVKAYVWDNNKD LAEW  
 LEKQLTEEDGVHSVIEENIKCISR DYVLKQIRSLVQANPEVAMDSI IHMTQHISPTQRAEVIRILS  
 TMDSPST **DYKDDDDKHHHHHH**

Black: ACC1  
 Red: FLAG-His-tags

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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)