



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Acc1

(Recombinant human protein with C-terminal His tag)

Catalog # **E-D017H**

Size: 10 µg

Product Description

Human ACC1 (acetyl-Coenzyme A carboxylase alpha transcript variant 1, GenBank Accession No. NM_198834), amino acids 39 – end with C-terminal His tag, MW = 266.5 kDa, expressed in baculovirus infected Sf9 cell expression system.

Purity

>20% by SDS-PAGE

Formulation

Recombinant protein is formulated in 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.1 mM EGTA, 0.1 mM EDTA, 10% glycerol and 3 mM DTT.

Storage and Stability

Store at or below -20° C. Enzyme will be stable for at least 6 months at -20° C as undiluted stock. Freeze in working aliquots to avoid repeated thawing and freezing.

Application

Acc1 is useful for the study of enzyme kinetics, regulation and dephosphorylation of target substrates.

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)