

Data Sheet

ACC, FLAG-Tag (*Candida albicans*)

Candida albicans, Recombinant, C-terminal
FLAG-tag

Catalog #: 100104

Lot #: 180817-A **Conc.:** 0.62 mg/ml

Description: *Candida albicans* ACC, also known as Acetyl-CoA carboxylase, GenBank Accession No. XM_713531, amino acids 1-2271 (end) with C-terminal FLAG-tag, expressed in Baculovirus infected Sf9 cell expression system. MW = 255 kDa.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, and 20% Glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Abu-Elheiga, L. *et al.*, *Science*. **291(5513)**, 2613-2616 (2001)
2. Tomas, E. *et al.*, *PNAS*. **99(25)**, 16309-16313. (2002)

Quality Assurance

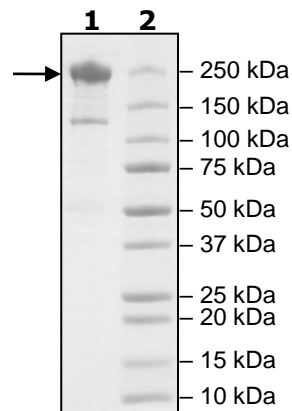
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg ACC

Lane 2:
Protein Marker

MW: 255 kDa

Purity: $\geq 79\%$



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