

Data Sheet

CEACAM, Avi-His-Tag, Biotin Labeled

Human, Recombinant, C-terminal Avi-His-Tag,
 Biotin Labeled

Catalog #: 70201

Lot #: 161020

Conc.: 0.93 mg/ml

Formulated in: 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C.

References:

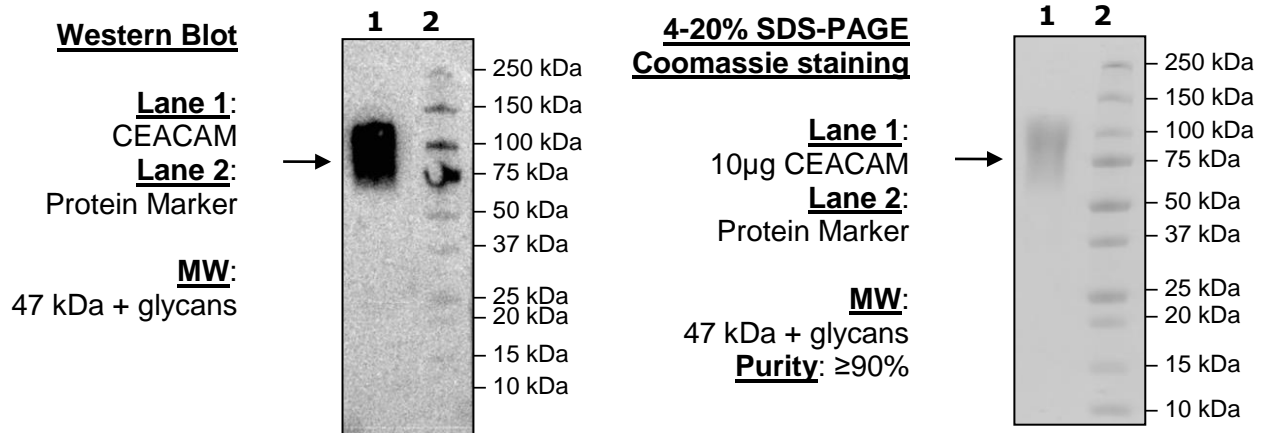
1. Carvalho, F. *et al.*, *J.E.M.* **206 (10)**, 2179 (2009).
2. Hauck R. *et al.*, *FEBS.* **478 (3)**, 260-266 (2000).

Description:

Human CEACAM, also known as Carcinoembryonic antigen-related cell adhesion molecule 1 or biliary glycoprotein, (GenBank Accession No. NM_001712), amino-acids 35-428, with C-terminal Avi-His-tag, MW = 47 kDa, expressed in HEK293 cells. This protein is enzymatically biotinylated using Avi-tag™ technology. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Application: Useful for western blot analysis.

Quality Assurance



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