

Data Sheet

CEACAM5, Avi-His-Tag

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

Catalog #: 100509

Lot#: 191209-1

Conc.: 1.49 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.

References:

1. Blumenthal, R.D., *et al.*, *BMC Cancer*. 2007; **7:2**
2. Saeland, E. *et al.*, *J. Cancer*. 2011; **131(1): 117-128**.

Description: Human CEACAM5, also known as Carcinoembryonic antigen-related cell adhesion molecule 5, CEA, and CD66e, GenBank Accession No. NM_004363.6, a.a. 35-685 with a C-terminal Avi-His-TagTM, expressed in a HEK293 cell expression system. MW=75 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Quality Assurance

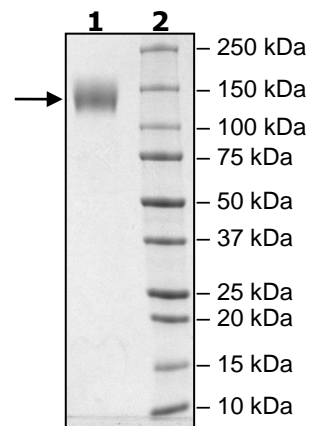
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg CEACAM5

Lane 2:
Protein Marker

MW: 75 kDa +glycans

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com