

Data Sheet

CRTAM, Fc-Fusion (IgG1), Avi-Tag

Human, Recombinant, C terminal Fc-Avi-Tag
Catalog #: 40686
Lot #: 160726-1 **Conc.:** 1.14 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 .*

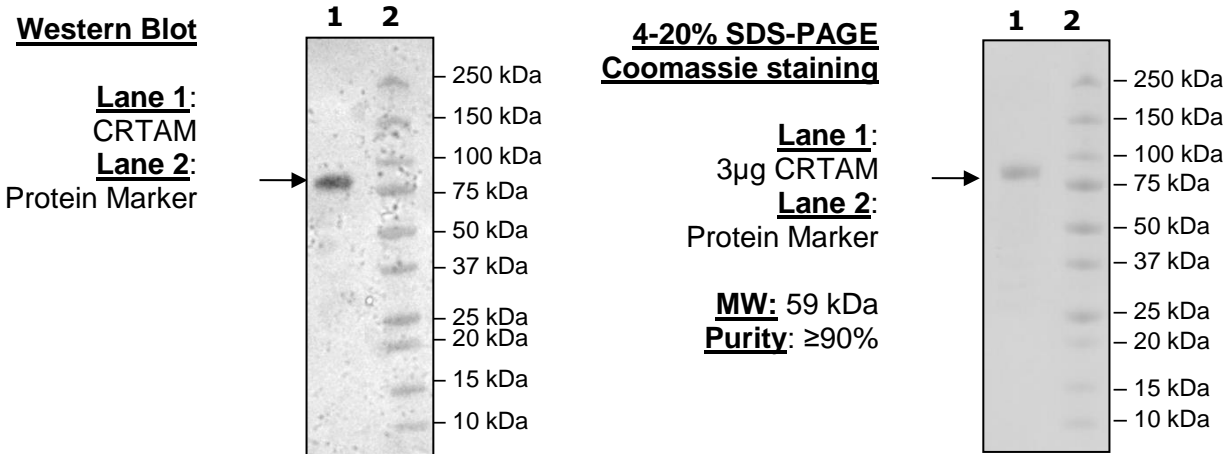
References:

1. Boles, K., *et al.*, *Blood*. 2005 Aug; **106(3)**: 779-86.
2. Yeh, J., *et al.*, *Cell*. 2008 Jan; **132(5)**:846-859.

Description: Human CRTAM, also known as Cytotoxic and Regulatory T-Cell Molecule, GenBank Accession No. NM_019604, amino acids 18-286, expressed in a HEK293 expression system. This protein has a C-terminal Avi-Tag fused to the Fc region of Human IgG1. MW = 59 kDa.

Application: Useful as a positive control for western blot studies.

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