

Data Sheet

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

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

Catalog #: 40687

Lot #: 170405

Conc.: 0.66 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.

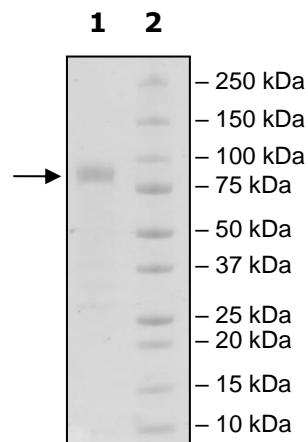
Description: Human CRTAM, also known as Cytotoxic and Regulatory T-Cell Molecule, GenBank Accession No. NM_019604, amino acids 18-286, expressed in a HEK293 expression system. This protein has a C-terminal Avi-Tag fused to the Fc region of Human IgG1 and is enzymatically biotinylated using Avi-tag™ technology.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

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg CRTAM
Lane 2:
Protein Marker

MW: 59 kDa
Purity: $\geq 86\%$



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