

Data Sheet

LAG3 (CD223)

Human, recombinant, Fc fusion protein
Catalog #: 71146
Lot#: 140422 Conc.: 1.43 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

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

Endotoxin Level: <0.1 EU/μg

References:

1. Byrnes, *et al.*. *Cytometry A*. 2012 Sep;**81(2)**:165-8.
2. Di, Carlo., *et al.*, *J. Pathol.* 2005 Jan; **205(1)**:82-91.

Description:

Human secreted LAG3, Fc fusion protein, also known as Lymphocyte-Activation Gene 3 and CD223. GenBank Accession No. NM_002286, a.a. 23-450 expressed in a HEK293 cell expression system. MW = 73.1 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful for studying protein binding and screening small molecules.

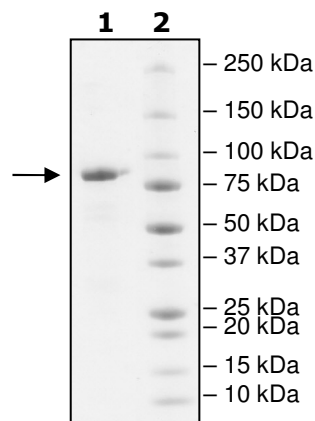
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg LAG3

Lane 2:
Protein Marker

MW: 73.1 kDa + glycans
Purity: ≥90%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

LQPGA EVPVWVAQEGAPAQLPCSPTIPLQDLSLLRRAGVTWQHQPDSGPPAAAPGH
PLAPGPHPAAPSSWGPRPRRYTVLSVGGGLRSGRLPLQPRVQLDERGRQRGDFSL
WLRPARRADAGEYRAAVHLRDRALSCRLRLRLGQASMTASPPGSLRASDWVILNCS
FSRPDRPASVHWFRNRGQGRVPVRESPPHHHLAESFLFLPQVSPMDSGPWGCILTYR
DGFNVSIMYNLTVLGLPPTPLTVYAGAGSRVGLPCRLPAGVGTSTRFLTAKWTPPGG
GPDLLVTGDNGDFTLRLEDVSAQAQAGTYTCHIHLEQQLNATVTLAIITVTPKSFGSP
GSLGKLLCEVTPVSGQERFVWSSLDTPSQRSFSGPWLEAQEAQLLSQPWQCQLYQ
GERLLGAAVYFTELSSPGAQRSGRAPGALPAGHLIEIEGRMDPKSCDKTHTCPPCPA
PELLGGPSVFLFPPKPKDTLMISRTPEVTCVVDVSHEDPEVKFNWYVDGVEVHNAKT
KPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCKVSNKALPAPIEKTISKAKGQPREPQ
VYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFL
YSKLTVDKSRWQQGNVFCSSVMHEALHNHYTQKSLSLSPGK

Black: LAG3 (a.a. 23-450)

Blue: linker

Red: Fc Fusion

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