

KAAG1 (human), (recombinant) (Fc-tag)

Ordering Information

[Order Online »](#)

ENZ-PRT330-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

kDa

M

116

66.2

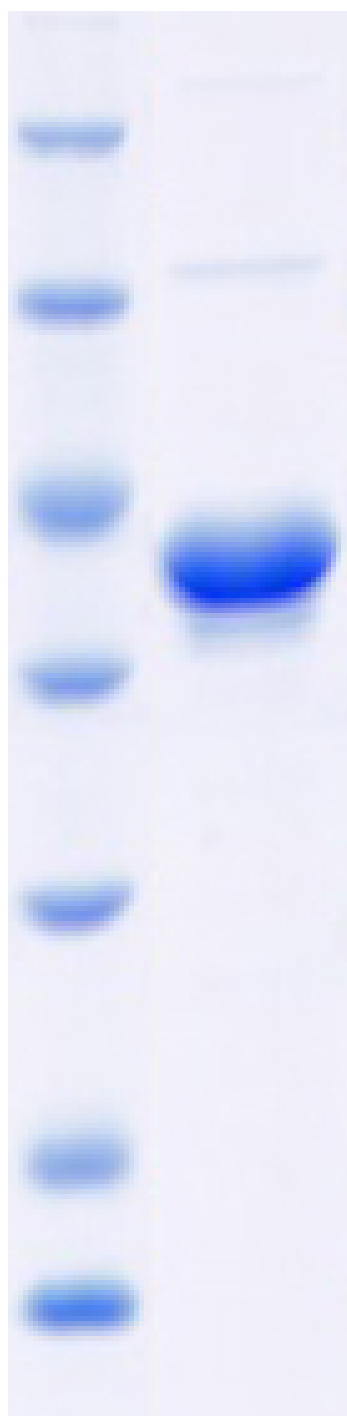
45.0

35.0

25.0

18.4

14.4



It migrates as approximately 25.1 kDa band in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Kidney-associated antigen 1, RU2AS, RU2 antisense gene protein

Endotoxin Content <1 EU/µg (LAL method)

Formulation Lyophilized from sterile 100 mM Glycine, 300 mM NaCl, 10% Glycerol, pH 8.2.

MW ~40.5 kDa (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in insect cells. Produced in a baculovirus expression system. A DNA sequence encoding the human KAAG1 (Met1-Lys84) was expressed with the Fc region of human IgG1 at the C-terminus.

UniProt ID Q9UBP8

Last modified: November 7, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)