

PRODUCT INFORMATION



Rabankyrin-5 (human, recombinant)

Item No. 45169

Overview and Properties

Synonyms: Ankyrin Repeat and FYVE Domain-containing Protein 1, Ankyrin Repeats Hooked To A Zinc Finger Motif, Rank-5

Source: Active recombinant human Rabankyrin-5 expressed in insect cells

Amino Acids: 2-1,169 (full length)

Uniprot No.: Q9P2R3

Molecular Weight: 128.4 kDa

Storage: -80°C (as supplied)

Stability: ≥1 year

Purity: ≥90% estimated by SDS-PAGE

Supplied in: 50 mM Tris-HCl, pH 7.5, with 200 mM sodium chloride and 20% glycerol

Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

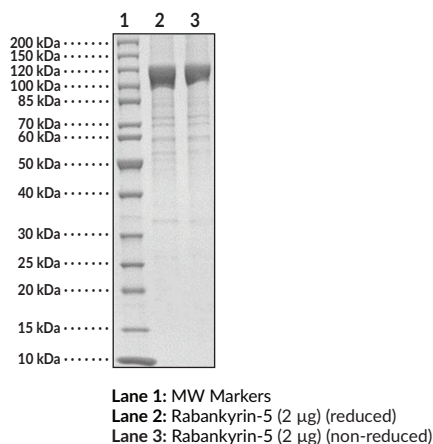
Protein Concentration: *batch specific* mg/ml

Activity: *batch specific* U/ml

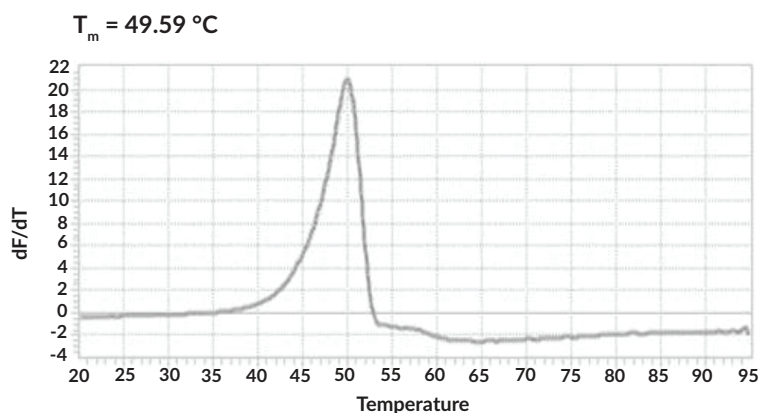
Specific Activity: *batch specific* U/mg

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images

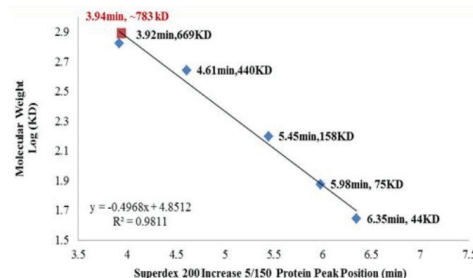
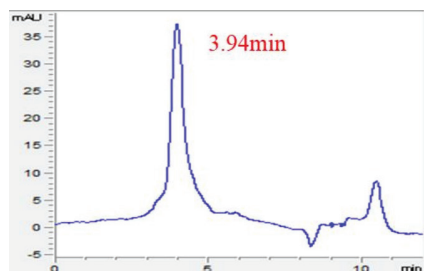


SDS-PAGE Analysis of Rabankyrin-5.



Protein melting thermal profile of Rabankyrin-5.

Analytical SEC



Analytical size exclusion chromatography of purified protein Rabankyrin-5.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

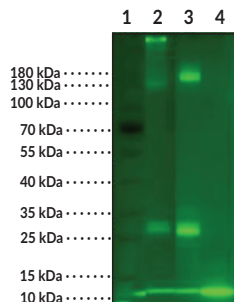
SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/09/2026

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Lane 1: MW Markers

Lane 2: BODIPY-Ub + UBE1 + UbcH5b + Rabankyrin-5 with ATP

Lane 3: BODIPY-Ub + UBE1 + UbcH5b with ATP

Lane 4: BODIPY-Ub + UBE1 + UbcH5b without ATP

Gel based ligase assay. The ubiquitin conjugating activity of Rabankyrin-5 was validated through its ability to catalyze the generation of polyubiquitin chains in the presence of the E1 activating enzyme UBE1, the E2 conjugating enzyme UbcH5b, and BODIPY-ubiquitin. Incubation of Rabankyrin-5 for 60 minutes at 37°C in the presence of BODIPY-ubiquitin, UBE1, UbcH5b and ATP (Lane 2) was compared alongside two control reactions with either Rabankyrin-5 (Lane 3) or Rabankyrin-5 with ATP (Lane 4) excluded from the reaction. Ubiquitin conjugates were identified by the migration of the Bodipy-ubiquitin band and these were observed only in the presence of ATP and Rabankyrin-5.

Description

Rabankyrin-5 is a Rab5 effector protein with roles in endosomal fusion, micropinocytosis, and mitochondrial homeostasis and dynamics.^{1,2} It is composed of an N-terminal BTB domain, 21 ankyrin (ANK) repeats, and a C-terminal FYVE finger domain and primarily localizes to endosomal membranes, but can also localize to the cytosol.¹ Rabankyrin-5 localizes with Rab5-positive early endosomes, and immunodepletion of Rabankyrin-5 using anti-Rabankyrin-5 antibodies inhibits early endosome fusion in HeLa cell cytosol. Knockdown of *ANKFY1*, the gene encoding Rabankyrin-5, using endoribonuclease-prepared small interfering RNA (esiRNA) inhibits EGF-induced micropinocytosis in skin cancer cells. *ANKFY1* knockdown also induces mitochondrial elongation and stasis in retinal pigment epithelial (RPE) cells.³ Cayman's Rabankyrin-5 (human, recombinant) protein has a calculated molecular weight of 128.4 kDa.

References

1. Schnatwinkel, C., Christoforidis, S., Lindsay, M.R., *et al.* The Rab5 effector Rabankyrin-5 regulates and coordinates different endocytic mechanisms. *PLoS Biol.* **2(9)**, E261 (2004).
2. Farmer, T., Naslavsky, N., and Caplan, S. Tying trafficking to fusion and fission at the mighty mitochondria. *Traffic* **19(8)**, 569-577 (2018).
3. Farmer, T., Reinecke, J.B., Xie, S., *et al.* Control of mitochondrial homeostasis by endocytic regulatory proteins. *J. Cell. Sci.* **130(14)**, 2359-2370 (2017).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM