

Product Information

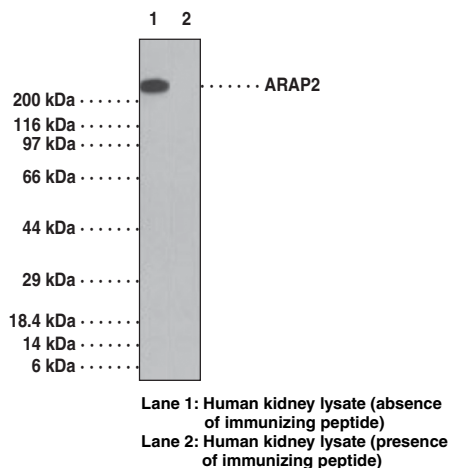


ARAP2 Polyclonal Antibody

Item No. 13495

- Contents:** This vial contains 100 µg of protein A-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Synonym:** Centaurin-δ-1
- Antigen:** Synthetic peptide from human ARAP2 within the region of amino acids 1,670-1,720
- Host:** Rabbit
- Cross Reactivity:** (+) Human ARAP2
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 1-2 µg/ml and 5 µg/ml for immunohistochemistry.

ARAP2, also known as centaurin-δ-1, is a protein with ARF-GAP, RHO-GAP, ankyrin repeat, RAS-associating, and pleckstrin homology domains. A phosphatidylinositol-3,4,5-trisphosphate (PtdIns-(3,4,5)-P₃)-dependent GTPase-activating protein, ARAP2, binds to RhoA-GTP, but lacks the predicted catalytic arginine in the RHO-GAP domain and does not have RHO-GAP activity. It modulates actin cytoskeleton remodeling by regulating ARF and RHO family members. Usually ARAP2 is activated by PtdIns-(3,4,5)-P₃ binding and can be activated by PtdIns-(3,4)-P₂ binding, although with lower efficiency. ARAP2 associates with focal adhesions and functions downstream of RhoA to regulate focal adhesion dynamics. It is usually detected in brain, thymus, lymph node, thyroid, spinal cord, trachea, heart, skeletal muscle, spleen, kidney, liver, placenta, lung, and peripheral blood leukocytes.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13495

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 04/12/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com