

Product Information

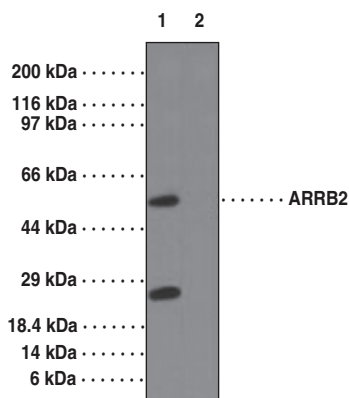


ARRB2 Peptide Affinity-purified Polyclonal Antibody

Catalog No. 13498

Contents:	This vial contains 100 µg of peptide affinity-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym:	β-Arrestin-2
Antigen:	Synthetic peptide from human ARRB2 within the region of amino acids 15-50
Host:	Rabbit
Cross Reactivity:	(+) Ovine, canine, equine, human, monkey, murine, and rat ARRB2
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	The recommended starting dilution for western blot is 0.25-0.5 µg/ml and 5 µg/ml for immunohistochemistry (paraffin-embedded sections).

β-Arrestin-2 (ARRB2) is an adaptor protein belonging to the arrestin family with an N-arrestin domain and a C-arrestin domain. ARRB2 is a regulatory protein involved in heterotrimeric G protein-coupled receptor desensitization and is known to regulate β-adrenergic receptor. A function, thus enhancing β2AR receptor mediated nuclear translocation of ERK. It binds to phosphorylated β-adrenergic receptors, thereby causing a significant impairment of their capacity to activate Gs proteins. Along with AIP4, ARRB2 acts as an endosomal sorting molecule that mediates CXCR4 entry into a degradative pathway. It may also be involved in hormone-specific desensitization of TSH receptors. ARRB2 is a Ca²⁺ binding protein of the retinal outer segments and binds to P-rhodopsin and is also involved in synaptic transmission in photoreceptor cells. It is widely expressed in most of the tissues, especially neuronal tissues and spleen.



Lane 1: HUVEC cell lysate (Absence of immunizing peptide)
Lane 2: HUVEC cell lysate (Presence of immunizing peptide)

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

MATERIAL 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 Material 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, 09/29/2009

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