

### Arf1 Activation Assay Kit ab211170

#### 2 图像

#### 概述

产品名称	Arf1 Activation Assay试剂盒
样品类型	Tissue, Adherent cells, Suspension cells
种属反应性	与反应: Mouse, Rat, Human
产品概述	Arf1 Activation Assay Kit (ab211170) uses GGA3 PBD Agarose beads to selectively isolate and pull-down the active form of Arf from purified samples or endogenous lysates. Subsequently, the precipitated GTP-Arf is detected by western blot analysis using an anti-Arf1 antibody, which detects the human, mouse and rat protein.

Features: 1) non radioactive assay format; 2) fast results: 1 hour assay plus electrophoresis/blotting time; 3) includes Cdc42 positive control; 4) pink colored agarose beads for easy identification during washing and aspiration steps.

#### 说明

Small GTP-binding proteins (or GTPases) are a family of proteins that serve as molecular regulators in signaling transduction pathways.

Arf1 (ADP-ribosylation factor 1) is a 21 kDa protein, member of ARF gene family, which constitutes part of the Ras superfamily. The Arf proteins are categorized as class I (Arf1, Arf2, Arf3), class II (Arf4 and Arf5) and class III (Arf6), and they all share a common gene organization. Arf1 regulates a variety of biological response pathways that include vesicle and protein trafficking, organelle structure, and endocytosis. Like other small GTPases, Arf1 cycles between an inactive GDP-bound form and an active GTP-bound form: in its active (GTP-bound) state, Arf1 binds specifically to the protein-binding domain (PBD) of GGA3 to control downstream signaling cascades.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

#### 性能

#### 存放说明

Store at -20°C. Please refer to protocols.

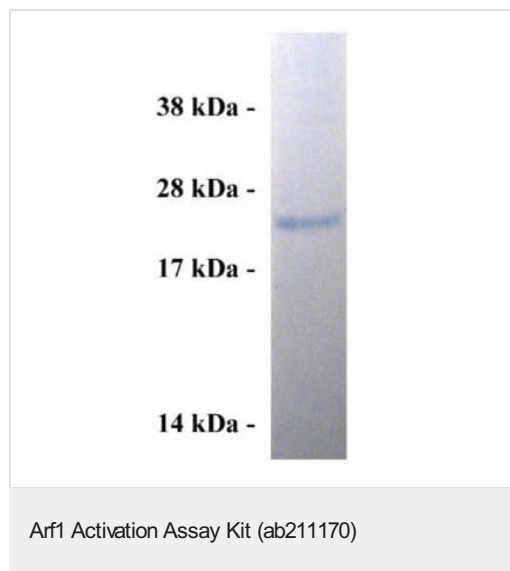
组件	20 tests
100X GDP	1 x 50µl
100X GTPγS	1 x 50µl
5X Assay/Lysis Buffer	1 x 30ml
ab211180 - GGA3 PBD Agarose Beads	1 x 800µl
anti-Arf1, Mouse monoclonal antibody	1 x 40µl
GTPase immunoblot Positive Control	1 x 100µl

功能	GTP-binding protein that functions as an allosteric activator of the cholera toxin catalytic subunit, an ADP-ribosyltransferase. Involved in protein trafficking among different compartments. Modulates vesicle budding and uncoating within the Golgi complex. Deactivation induces the redistribution of the entire Golgi complex to the endoplasmic reticulum, suggesting a crucial role in protein trafficking. In its GTP-bound form, its triggers the association with coat proteins with the Golgi membrane. The hydrolysis of ARF1-bound GTP, which is mediated by ARFGAPs proteins, is required for dissociation of coat proteins from Golgi membranes and vesicles.
序列相似性	Belongs to the small GTPase superfamily. Arf family.
细胞定位	Golgi apparatus. Cytoplasm > perinuclear region.

图片



Arf1 activation assay. Lane 1: HeLa cell lysate loaded with GDP and incubated with GGA3 PBD Agarose beads. Lane 2: HeLa cell lysate loaded with GTPγS and incubated with GGA3 PBD Agarose beads.



Arf1 positive control (293 cell lysate at 0.5 mg/mL) detected by immunoblot.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

#### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

#### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors