V. 1.0

Arp2/3 Protein Complex Source: Bovine brain Cat. # RP01

Upon arrival store at 4°C (desiccated)
See datasheet for storage after reconstitution

Material

The Arp2/3 complex (Actin-Related Proteins) has been purified from bovine brain and consists of seven evolutionarily conserved protein subunits (1), all present in approximately equal stoichiometry (see Figure 1). The Arp2/3 complex (molecular weight of 224 kDa) is a key regulator of branched actin filament nucleation and is important for cell locomation, phagocytosis and intracellular motility. Arp2/3 complex is supplied as a white lyophilized powder.

Storage and Reconstitution

Store lyophilized product desiccated (<10% humidity) at 4°C where it is stable for 6 months. Resuspend the protein complex to 5 mg/ml with 10 μ l of cold Milli-Q water. When resuspended the complex is in the following buffer: 20 mM Tris pH 7.5, 25 mM KCl, 1 mM MgCl $_2$, 0.5 mM EDTA, 0.1 mM ATP, 1.0% (v/v) dextran and 5% (v/v) sucrose. The protein should then be aliquoted into experiment sized amounts, snap frozen in liquid nitrogen and stored at -70°C where it is stable for 6 months. Avoid multiple freeze-thaw cycles. Further dilution of Arp2/3 should be made in the following buffer: 20 mM Tris pH 7.5, 25 mM KCl, 1 mM MgCl $_2$ and 1 mM DTT (Note: add DTT to the buffer immediately prior to use).

Purity

Protein purity is determined by scanning densitometry of Coomassie Blue stained protein on a 4-20% polyacrylamide gel. Arp2/3 protein complex is determined to be 90% pure (see Figure 1).

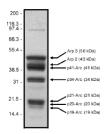


Figure 1. Arp2/3 Protein Purity Determination. A 10 μg sample of Arp2/3 complex was separated by electrophoresis in a 4-20% SDS-PAGE system and stained with Coomassie Blue. Protein quantitation was determined with the Precision RedTM Protein Assay Reagent (Cat. # ADV02). Mark12 molecular weight markers are from Invitrogen.

Biological Activity Assay

The Arp2/3 complex is able to induce the branched polymerization of actin filaments *in vitro* at a molar ratio of 1:200 (Arp2/3:actin). This stimulation is observable in an *in vitro* polymerization assay (described below), however, the stimulation from Arp2/3 alone is very low under these conditions. In the presence of N-WASP protein (or the VCA domain of N-WASP, Cat. # VCG03) the nucleating activity of Arp2/3 is greatly enhanced. The *in vitro* polymerization assay is described below:

Reagents

- 1. Arp2/3 protein complex (Cat. # RP01)
- VCA domain-GST fusion (Cat # VCG03)
- 3. Pyrene labeled actin (Cat. # AP05)
- Polymerization buffer 1.5x stock: 7.5 mM Tris pH 7.5, 75 mM KCl, 3 mM MgCl₂, 1.5 mM EGTA, 0.15 mM CaCl, 0.75 mM DTT, 0.3 mM ATP [add fresh from a 100 mM stock pH 7.0, immediately prior to use].
- General Actin Buffer (Cat. # BSA01):10 mM Tris pH 7.5, 0.2 mM CaCl₂, 0.2 mM ATP, 1 mM DTT.

Equipment

- Fluorescence spectrophotometer with an excitation wavelength of 360 +/- 20 nm and an emission wavelength of 405 +/- 10 nm or 420 +/- 20 nm. Cytoskeleton recommends the SPECTRAFluor Plus from TECAN Austria GmbH or Gemini from Molecular Devices Inc.
- Black polystrene 96 well assay plate (Costar, Cat. # 3915).

Method

- Resuspend and dilute pyrene labeled muscle actin (Cat. # AP05) to 0.45 mg/ml (10.4 µM) with General Actin Buffer (Cat. # BSA01) supplemented with 0.2 mM ATP and 1 mM DTT. Leave on ice for 1 h to depolymerize actin oligomers.
- Centrifuge the actin at 14,000 rpm at 4°C for 30 min to remove residual nucleating centers.
- Pipette the top 80% of the supernatant into a new microfuge tube on ice.
- Dilute the Arp2/3 complex (Cat. # RP01) to 0.3 mg/ml in Gbuffer. Keep on ice.
- Resuspend one tube of VCA domain protein (Cat. # VCG03) to 1 mg/ml by adding 500 µl of Milli-Q water. Pipet up and down slowly to resuspend the pellet. Keep on ice.
- Just before use, dilute the pyrene-labeled actin to 0.1 mg/ ml in ice-cold General Actin Buffer.

7. Add the following components to the 96 well assay plate:

Well	Poly buffer (μl)	RP01 (µI)	VCA domain (μl)
A1	200	0	0
B1	200	0	0
C1	200	0	5
D1	200	0	5
E1	200	2	0
F1	200	2	0
G1	200	2	5
H1	200	2	5

- Using a multi-channel pipet, add 100 µl of diluted pyrenelabeled actin to wells A1-H1 of the assay plate. Note: Do not introduce air bubbles into the wells.
- Place the 96 well plate into the fluorescent spectrophotometer and read the samples for 1 h.
- In the assay described above, pyrene actin is present at a final concentration of 0.8 µM, Arp2/3 complex at 10 nM and VCA domain at 400 nM.
- Results for a typical actin polymerization assay is shown in Figure 2.

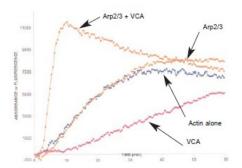


Figure 2. Actin Polymerization Assay with Arp2/3 and VCA domain proteins. Actin polymerization was carried out as described in the method; all reactions contain pyrene labeled actin. Reactions containing Arp2/3 or VCA in addition to actin show little or no enhancement of actin nucleation. In the presence of the VCA domain however, Arp2/3 results in an enhancement of actin nucleation (Note the steep nucleation phase of polymerization). Actin polymerization is measured in arbitrary fluorescent units over time.

Product Uses

- Characterization of actin polymerization factors
- · Formation of branched actin filaments
- Antibody standard for Western blot analysis

References

 Pollard, T.D. et al. 2000. Ann. Rev. Biophys. Biomol. Struct. 29: 545-576.

Product Citations/Related Products

For the latest citations and related products please visit www.cytoskeleton.com.