

Recombinant ACF complex

Catalog No: 31509, 31913

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: *Drosophila melanogaster*

Buffer Contents: Recombinant ACF complex is supplied in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP. Please refer to product insert upon arrival for lot-specific concentration.

Background: ACF Complex (ATP-utilizing chromatin assembly and remodeling factor protein complex) which contains Acf1 and Iswi catalyzes the ATP-dependent assembly of periodic nucleosome arrays in vitro. The catalytic subunit of ACF is the ISWI ATP-dependant chromatin remodeling protein. ACF is related to CHRAC (Chromatin Accessibility Complex), which also contains Acf1 and Iswi. ACF is involved in the establishment and possibly maintenance of silenced heterochromatic domains in the genome, both at pericentric heterochromatin and at genes repressed by the Polycomb complex.

Protein Details: Recombinant ACF complex that includes full length *Drosophila* Acf1 and Iswi (accession numbers NP_536734.2 and NP_523719.1, respectively) was expressed in Sf9 cells and contains an FLAG-Tag at the N-terminus of Acf1. The molecular weights of expressed Acf1 and Iswi are 172 kDa and 119 kDa, respectively.

Application Notes: Recombinant ACF complex is suitable for in vitro chromatin assembly. This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

References:

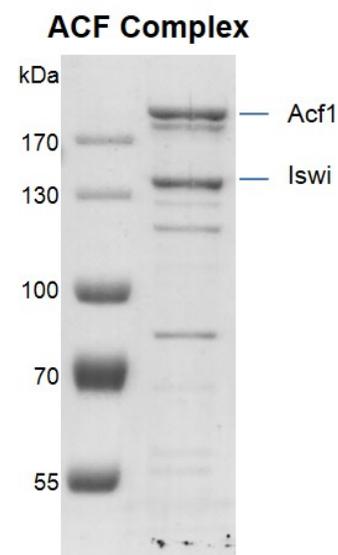
This product was used in the following publications:

Mol. Cell (2018). 69(3): 385-397. PMID: 29336876.

Cell (2019). 177(7): 1781-1796. PMID: 31104845.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant ACF complex gel

ACF complex was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.

MW: Acf1: 172 kDa. Iswi: 119 kDa.

Purity: > 75%.