

ACF Chromatin Remodeling Enzyme Complex



EpiCypher®

Catalog No 15-1013
Lot No 18200001
Pack Size 100 reactions

Type Remodeler/ATPase **Expressed In** SF9 cells
Mol. Wgt. 185 kDa **Epitope Tag** 6His-FLAG

Product Description:

Recombinant *Drosophila* ACF Chromatin Remodeling Enzyme Complex, assembled from Acf1-FLAG and ISWI produced in SF9 cells. ACF is an ATP-dependent chromatin remodeling complex that regulates nucleosome spacing.

Formulation:

ACF Complex at 0.18 mg/mL in 5.2 μ L of 20 mM Tris HCl pH 7.9, 150 mM NaCl, 2 mM MgCl₂, 0.2 mM EDTA, 10 mM beta-glycerophosphate, 0.01% NP-40, 0.2 mM PMSF, 0.5 mM benzamidine, 1 mM DTT, 15% glycerol, 2 μ g/mL leupeptin, 1 μ g/mL aprotinin, 0.4 mg/mL recombinant insulin, 0.4 mg/mL 3x FLAG peptide. Molarity = 973 nM

Storage and Stability:

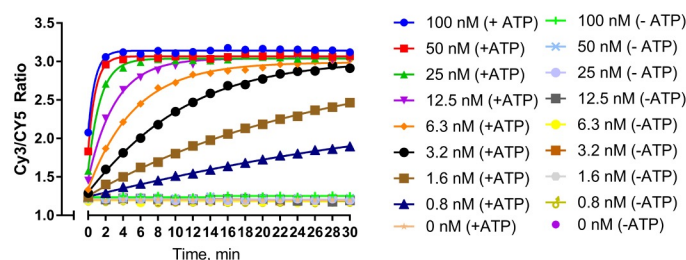
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

This product is sufficient to perform 100 remodeling reactions using EpiDyne®-FRET substrate. A single reaction is defined as 10 μ L containing [5 nM ACF Complex, 20 nM EpiDyne®-FRET (Catalog No. 16-4201), 50 μ M rATP] and remodels to completion in <20 minutes. 5x ACF Remodeling Assay Buffer is included (Catalog No. 21-0013; 100 mM Tris HCl pH 7.5, 250 mM KCl, 15 mM MgCl₂, 0.05% (w/v) BSA, 0.05% (v/v) Tween 20). For more information, see the EpiDyne®-FRET Technote (<https://www.epicypher.com/resources/technical-notes/>) or contact techsupport@epicypher.com

References:

Fyodorov DV & Kadonaga JT (2002) *Nature* 418: 897-900.



ATP-dependent Chromatin Remodeling Assay: EpiDyne®-FRET Chromatin Remodeling Substrate (20 nM; Catalog No. 16-4201) incubated with ACF Remodeling Enzyme Complex (concentrations indicated), with or without ATP (50 μ M) in 1x ACF Remodeling Assay Buffer. Curves denote FRET efficiency/chromatin remodeling.

This product is for *in vitro* research use only and is not intended for use in humans or animals.

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com