

# 5x ACF Remodeling Assay Buffer



EpiCypher®

**Catalog No** 21-0013  
**Lot No** 20288002-39  
**Pack Size** 500  $\mu$ L

## Product Description:

5x concentrated reaction buffer for use with ACF Remodeling Enzyme Complex (EpiCypher 15-1013). Optimized for use with recombinant *Drosophila* ACF 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:

5x ACF Remodeling Assay Buffer (100 mM Tris HCl pH 7.5, 250 mM KCl, 15 mM MgCl<sub>2</sub>, 0.05% (w/v) BSA, 0.05% (v/v) Tween 20) is formulated for use with ACF Remodeling Enzyme Complex.

## Storage and Stability:

For best results, store 5x ACF Remodeling Assay buffer at 4°C and avoid freezing.

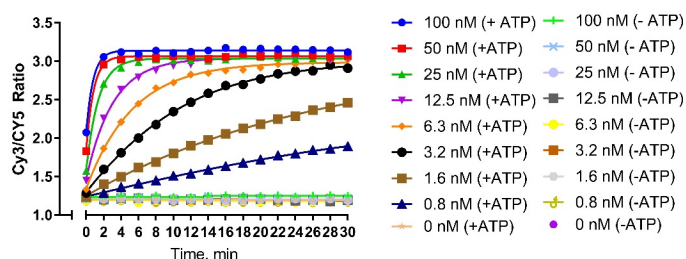
## Application Notes:

This product is formulated to perform remodeling reactions using ACF Remodeling Enzyme Complex and EpiDyne®-FRET substrate in a reaction defined as 10  $\mu$ L containing [5 nM ACF Complex, 20 nM EpiDyne-FRET, and 50  $\mu$ M rATP] and remodels to completion in <30 minutes.

For further information, please contact [techsupport@epicypher.com](mailto:techsupport@epicypher.com) or refer to the EpiDyne-FRET technote at [www.epicypher.com/technotes](http://www.epicypher.com/technotes).

## References:

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



**ATP-dependent Chromatin Remodeling Assay:** EpiDyne-FRET Chromatin Remodeling Substrate (20 nM; EpiCypher 16-4201) incubated with ACF Remodeling Enzyme Complex (EpiCypher 15-1013) using ACF Remodeling Assay Buffer with or without ATP (50  $\mu$ M). Curves denote FRET efficiency/chromatin remodeling.