

Nucleosome Remodeling Assay by EpiDyne™-FRET

Fully recombinant nucleosome substrates engineered for chromatin remodeling studies

Authors:

Matthew F. Whelihan, Martis W. Cowles,
Brandon A. Boone, Zu-Wen Sun, James R. Bone,
Michael-Christopher Keogh

EpiCypher Inc. Research Triangle Park, NC 27713
Correspondence: info@epicypher.com

Overview

Chromatin remodeling, or the repositioning of nucleosomes, regulates DNA access and thus gene expression and genome repair. Many ATP-dependent remodeling enzyme complexes are associated with human disease but are challenging study targets due to the requirement for nucleosome-based substrates. EpiCypher has addressed this need by developing the EpiDyne platform of fully recombinant remodeling substrates to monitor nucleosome repositioning along DNA. Here, we demonstrate the utility of EpiDyne-FRET in a homogeneous remodeling assay using Fluorescence Resonance Energy Transfer (FRET) readout.

Chromatin Remodeling Enzymes As Therapeutic Targets

Aberrant nucleosome organization can severely disrupt gene expression, DNA repair and cellular differentiation, and it also plays a major role in human disorders, including cancer, inflammation, autoimmunity, schizophrenia, cardiovascular disease, and intellectual disability. Remarkably, nearly 20% of all cancers contain mutations in subunits from the SWI/SNF family of ATP-dependent chromatin remodeling complexes. These enzyme complexes regulate local genome access by 'pumping' the DNA around histone octamers, thus 'sliding' nucleosomes [1]. Recurrent somatic mutations in SWI/SNF subunits are observed in multiple cancers, supporting a driver role in tumorigenesis [2]. The mutated remodeling proteins are attractive therapeutic targets, since further compromising their ATPase activity promotes cancer cell death but spares normal cells [2, 3]. This phenomenon is known as synthetic lethality, and identifying inhibitors to exploit it may lead to drugs with cancer specificity [4, 5].

Chromatin Remodeling Assays Using Recombinant Mononucleosome Substrates

EpiCypher has developed EpiDyne-FRET to study the function of chromatin remodeling enzyme complexes. The nucleosome consists of a core histone octamer wrapped in ~147 bp DNA and represents the basic repeating unit of chromatin. EpiDyne-FRET is comprised of a terminally positioned histone octamer (H2A-T120C*Cy5) wrapped in 5' Cy3-labelled DNA [8] (Figure 1). In its assembled start position, Cy3-Cy5 FRET is at a maximum, with nucleosome remodeling detected by loss of the Cy5 signal as the histone octamer is relocated (towards the template DNA 3' end) or ejected.

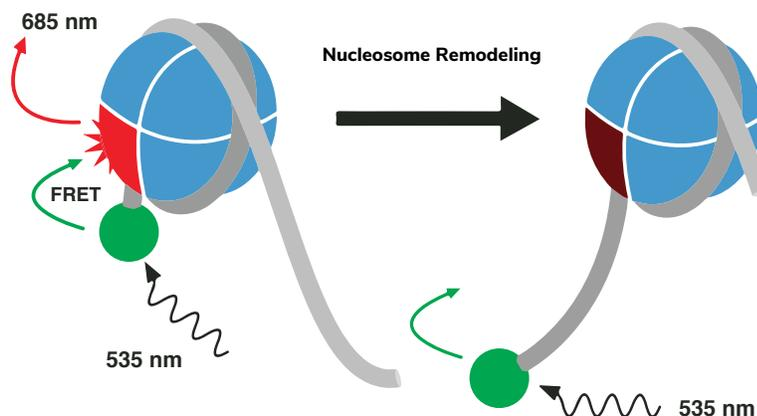


FIGURE 1: EPIDYNE-FRET NUCLEOSOME REMODELING SUBSTRATES consist of a Cy5-labeled human histone octamer (H2A T120C-Cy5; shown as red section of octamer) wrapped by 5' Cy3-labeled DNA (217bp; green ball) comprising a terminally nucleosome positioning sequence (147bp Widom 601 [7]) adjacent to a TGGA-repeat region refractory to nucleosome assembly [8]. In its assembled starting state, Cy3-Cy5 FRET is at a maximum. The activity of an ATP-dependent remodeler (e.g. RSC or another SWI/SNF ATPase) is detected by a reduction in FRET signal as the Cy3-labeled DNA 5' end is moved away from the Cy5-labeled octamer. EpiDyne-FRET is a one-step no-wash method immediately compatible with HTS applications.

Continued...

Nucleosome Remodeling Assay by EpiDyne™-FRET

Fully recombinant nucleosome substrates engineered for chromatin remodeling studies

Continued from page 1

Reagents and Materials Required for the Assay [see Note (i)]

- Yeast RSC [9] (substrate also compatible with dACF [ACF-ISWI], dNoRC [Toutatis-ISWI] and human SMARCA2/4 (BRG1/BRM); data not shown)
- EpiDyne-FRET nucleosome remodeling substrate (EpiCypher Catalog No. 16-4201)
- Assay Buffer (20 mM Tris pH 7.5, 50 mM KCl, 3 mM MgCl₂ and 0.1 mg/mL BSA (Sigma Catalog No. A3059))
- ATP (Invitrogen Catalog No. PV3227)
- Optional: ATPγS (a non-hydrolyzable form of ATP)
- Corning 3820 384-well assay plate (Fisher Catalog No. 07-200-891)
- 16 channel multipipetter for 384-well plates
- 384-well Fluorescence microplate reader capable of Cy3/Cy5-FRET detection (e.g. Envision, Perkin Elmer)

Standard Protocol

1. Determine the amount of RSC (10 nM final), EpiDyne-FRET (20 nM final), and ATP (2 mM final) needed - based on the intended time points / replicates (triplicate shown) [see Notes (ii & iii)].
2. Prepare reaction components (room temp).
 - i. 2x Enzyme/Substrate solution (20 nM RSC + 40 nM EpiDyne-FRET substrate in 2x Assay Buffer)
 - ii. 2x ATP +/- inhibitor solution(s) (4 mM ATP in ddH₂O + inhibitors [e.g. ATPγS] as required)
3. Add 5 μL of 2x Enzyme/Substrate solution to the microplate.
4. Initiate reactions by adding 5 μL of the 2x ATP +/- inhibitor solution.
5. At appropriate timepoints read in a 384-well plate reader capable of Cy3 (excitation-531 nm / emission-579 nm) / Cy5 (emission-685 nm) detection.
6. Data is expressed as the ratio of the raw Cy3 and Cy5 emission signals at each time point [8].

NOTES

- (i) Store protein aliquots at -80°C and avoid freeze/thaw.
- (ii) For assay development it is advised to titrate the concentration of multiple reaction components, including enzyme, substrate, and ATP.
- (iii) Since data is expressed as Cy3-Cy5 emission it is important not to add excess substrate to these reactions.
- (iv) Reactions can begin rapidly. If setting up a large plate without an auto-injector there is no capability of a T0 for each condition.

References

1. Hamiche, A., et al., ATP-dependent histone octamer sliding mediated by the chromatin remodeling complex NURF. *Cell*, 1999. 97(7): p. 833-42.
2. Vangamudi, B., et al., The SMARCA2/4 ATPase Domain Surpasses the Bromodomain as a Drug Target in SWI/SNF-Mutant Cancers: Insights from cDNA Rescue and PFI-3 Inhibitor Studies. *Cancer Res*, 2015. 75(18): p. 3865-78.
3. Wu, Q., et al., Targeting the chromatin remodeling enzyme BRG1 increases the efficacy of chemotherapy drugs in breast cancer cells. *Oncotarget*, 2016. 7(19): p. 27158-75.
4. Helming, K.C., X. Wang, and C.W. Roberts, Vulnerabilities of mutant SWI/SNF complexes in cancer. *Cancer Cell*, 2014. 26(3): p. 309-17.
5. Karnezis, A.N., et al., Dual loss of the SWI/SNF complex ATPases SMARCA4/BRG1 and SMARCA2/BRM is highly sensitive and specific for small cell carcinoma of the ovary, hypercalcaemic type. *J Pathol*, 2016. 238(3): p. 389-400.
6. Clapier, C.R., et al., Regulation of DNA Translocation Efficiency within the Chromatin Remodeler RSC/ Sth1 Potentiates Nucleosome Sliding and Ejection. *Mol Cell*, 2016. 62(3): p. 453-61.
7. Lowary, P.T. and J. Widom, New DNA sequence rules for high affinity binding to histone octamer and sequence-directed nucleosome positioning. *J Mol Biol*, 1998. 276(1): p. 19-42.
8. Cao, H., et al., TGGA repeats impair nucleosome formation. *J Mol Biol*, 1998. 281(2): p. 253-60.

Acknowledgments

We would like to thank Drs. Cedric Clapier and Bradley Cairns for kindly providing recombinant RSC complex used for this study. Development for this product was sponsored by National Institute Of General Medical Sciences of the National Institutes of Health (Award# 1R43GM123869-01)

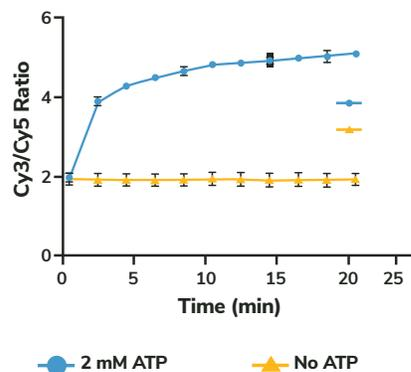


FIGURE 2: RSC/ATP-DEPENDENT NUCLEOSOME REMODELING REACTION. EpiDyne-FRET nucleosomes (20 nM) were incubated with RSC chromatin remodeler (10 nM) in the presence or absence of 2 mM ATP (conditions previously determined as optimal for this enzyme; data not shown). Upon ATP addition, reactions were read in an Envision Multi-label plate reader at times indicated. Data is presented as the mean±SD of the Cy3/Cy5 ratio (N=6) [see Note (iv)].

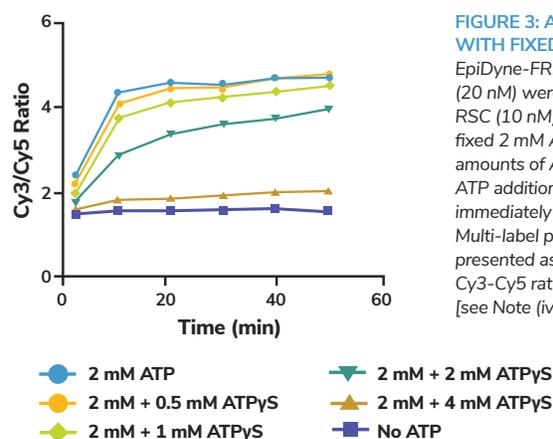


FIGURE 3: ATPγS TITRATION WITH FIXED ATP. EpiDyne-FRET nucleosomes (20 nM) were incubated with RSC (10 nM) in the presence of fixed 2 mM ATP with increasing amounts of ATPγS. Upon ATP addition, reactions were immediately read in an Envision Multi-label plate reader. Data is presented as the mean of the Cy3-Cy5 ratio (N=2) [see Note (iv)].

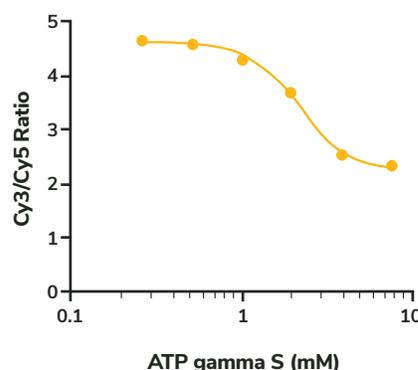


FIGURE 4: IC50 CALCULATION FOR ATPγS. Data from Figure 3 (20-minute time point) were analyzed to determine the IC50 value for ATPγS (2.29 mM). Data is presented as the mean of the Cy3-Cy5 ratio vs. the ATPγS concentration.