

# 5x SMARCA Remodeling Assay Buffer



## EpiCypher®

**Catalog No** 21-0014  
**Lot No** 20288002-16  
**Pack Size** 500  $\mu$ L

### Product Description:

5x concentrated reaction buffer for use with SMARCA Remodeling Enzymes (EpiCypher 15-1014 and 15-1015). Optimized for use with SMARCA Remodeling Enzyme, produced in SF9 cells. SMARCA is an ATP-dependent chromatin remodeling complex that regulates nucleosome spacing.

### Formulation:

5x SMARCA 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 SMARCA Remodeling Enzyme.

### Storage and Stability:

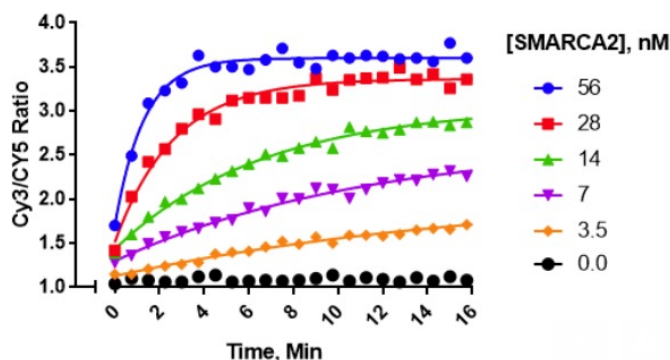
For best results, store 5x SMARCA Remodeling Assay buffer at 4°C and avoid freezing.

### Application Notes:

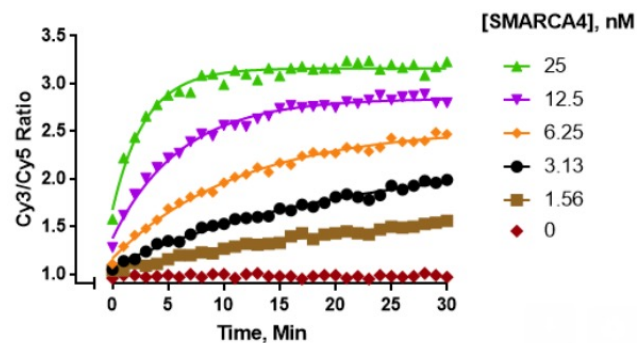
This product is formulated to perform remodeling reactions using SMARCA Remodeling Enzyme and EpiDyne®-FRET substrate (EpiCypher 16-4201) in a 10  $\mu$ L reaction containing 20 nM SMARCA, 20 nM EpiDyne-FRET, and 2.5 mM rATP. Enzyme remodels to completion in <30 minutes.

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:



**ATP-dependent Chromatin Remodeling Assay:** EpiDyne-FRET Chromatin Remodeling Substrate (20 nM; EpiCypher 16-4201) incubated with SMARCA2 Remodeling Enzyme (EpiCypher 15-1015; concentrations indicated). Curves denote FRET efficiency and chromatin remodeling.



**ATP-dependent Chromatin Remodeling Assay:** EpiDyne-FRET Chromatin Remodeling Substrate (20 nM; EpiCypher 16-4201) incubated with SMARCA4 Remodeling Enzyme (EpiCypher 15-1014; concentrations indicated). Curves denote FRET efficiency and chromatin remodeling.

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