

5x SMARCA Remodeling Assay Buffer



EpiCypher®

Catalog No 21-0014
Lot No 20288002-16
Pack Size 500 μ 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 MgCl₂, 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.

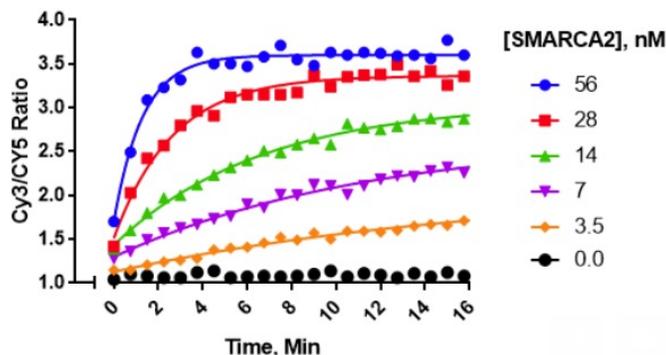
Application Notes:

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

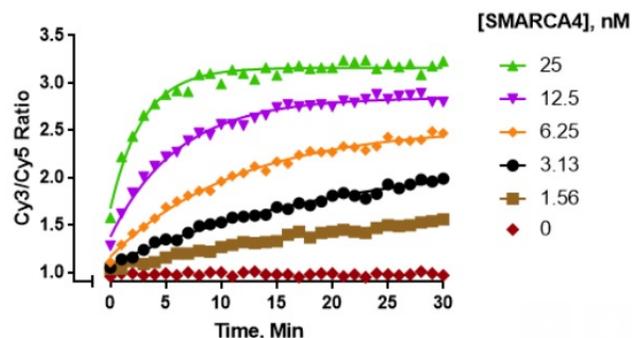
For further information, please contact techsupport@epicypher.com or refer to the EpiDyne-FRET technote at www.epicypher.com/technotes.

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 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.