

# SMARCA2 Chromatin Remodeling Enzyme



EpiCypher®

**Catalog No** 15-1015  
**Lot No** 22073006-01  
**Pack Size** 100 Reactions

**Type** Remodeler/ATPase      **Expressed In** SF9 cells

**Mol. Wgt.** 181 kDa      **Epitope Tag** FLAG

## Product Description:

Enzymatically active, full length recombinant human SMARCA2 Remodeling Enzyme (BRM, UniProt reference no. P51531-2) produced in SF9 cells. SMARCA2 is the catalytic subunit of the SWI/SNF complex. It is an ATP-dependent chromatin remodeling enzyme that regulates nucleosome spacing.

## Formulation:

SMARCA2 at 0.26 mg/mL in 13.9  $\mu$ L of 25 mM HEPES pH 7.6, 300 mM NaCl, 0.1 mM EDTA, 10% glycerol, 1 mM DTT, 0.4 mM PMSF, 0.1 mM benzamidine, 0.4 mg/mL recombinant human insulin. Molarity = 1.44  $\mu$ M.

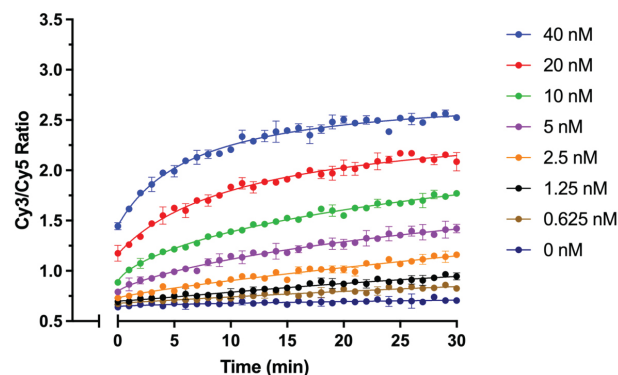
## Storage and Stability:

Stable for six months at  $-80^{\circ}\text{C}$  from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

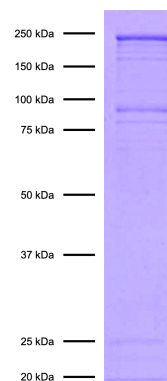
## Application Notes:

This product is sufficient to perform 100 remodeling reactions using EpiDyne®-FRET substrate (EpiCypher 16-4201). A single reaction is defined as 10  $\mu$ L containing [20 nM SMARCA2, 20 nM EpiDyne®-FRET, 1 mM rATP] and remodels to completion in <30 minutes. 5x SMARCA Remodeling Assay Buffer is included (EpiCypher 21-0014). For more information, see the EpiDyne®-FRET Technote ([epicypher.com/resources/technical-notes](http://epicypher.com/resources/technical-notes)) or contact [techsupport@epicypher.com](mailto:techsupport@epicypher.com).

## References:



**Figure 1: ATP-dependent chromatin remodeling assay.** EpiDyne®-FRET Chromatin Remodeling Substrate (EpiCypher 16-4201; 15nM) incubated with SMARCA2 Remodeling Enzyme (concentrations indicated) in 1x SMARCA2 Remodeling Assay Buffer. Curves denote FRET efficiency/chromatin remodeling.



**Figure 2: Protein gel data.** SMARCA2 Remodeling Enzyme (1  $\mu$ g) was run on an SDS-PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

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