

SMARCA4 Chromatin Remodeling Enzyme



EpiCypher®

Catalog No 15-1014
Lot No 21145007-02
Pack Size 100 reactions

Type Remodeler/ATPase **Expressed In** SF9 cells
Mol. Wgt. 186.6 kDa **Epitope Tag** 6xHis/FLAG

Product Description:

Full length recombinant human SMARCA4 Remodeling Enzyme (BRG1, UniProt reference no. P51532-1) produced in SF9 cells. SMARCA4 is a crucial component of the SWI/SNF complex. It is an ATP-dependent chromatin remodeling enzyme that regulates nucleosome spacing.

Formulation:

SMARCA4 at 0.13 mg/mL in 14.4 μ L of 25 mM HEPES pH 7.6, 0.1 mM EDTA, 10% glycerol, 300 mM NaCl, 1 mM DTT, 0.4 mM PMSF, 1 mM benzamidine, 0.4 mg/mL insulin. Molarity = 0.697 μ M

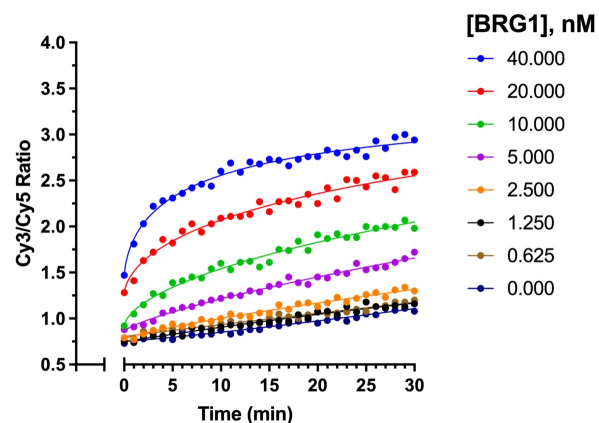
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

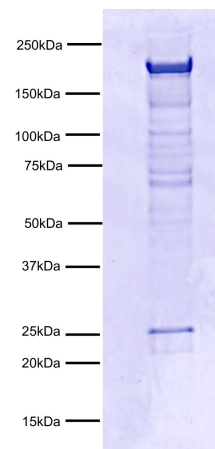
Application Notes:

This product is sufficient to perform 100 remodeling reactions using EpiDyne®-FRET substrate (Catalog No. 16-4201). A single reaction is defined as 10 μ L containing [10 nM SMARCA4, 20 nM EpiDyne®-FRET, 1 mM rATP] and remodels to completion in <30 minutes. 5x SMARCA Remodeling Assay Buffer is included (Catalog No. 21-0014; 100 mM Tris HCl pH 7.5, 250 mM KCl, 15 mM MgCl₂, 0.05% (w/v) BSA, 0.05% (v/v) Tween 20). For more information, see the EpiDyne®-FRET Technote (<https://www.epicypher.com/resources/technical-notes/>) or contact techsupport@epicypher.com

References:



ATP-dependent Chromatin Remodeling Assay: EpiDyne®-FRET Chromatin Remodeling Substrate (20 nM; Catalog No. 16-4201) incubated with SMARCA4 Remodeling Enzyme (concentrations indicated) in 1x SMARCA4 Remodeling Assay Buffer. Curves denote FRET efficiency/chromatin remodeling.



Protein Gel Data: SMARCA4 Remodeling Enzyme (1 μ 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.