

SMARCA4 Chromatin Remodeling Enzyme

Catalog No. 15-1014
Lot No. 18320001
Pack Size 100 reactions



EpiCypher®

Type Remodeler/ATPase **Expressed In** SF9 cells
Mol. Wgt. 182.7 kDa **Epitope Tag** FLAG

Product Description:

Full length recombinant human SMARCA4 (BRG1) Remodeling Enzyme 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 0.85 mg/ml in 4.3 µl of 20 mM Tris HCl pH 7.9, 150 mM NaCl, 2 mM MgCl₂, 0.2 mM EDTA, 10 mM beta-glycerophosphate, 0.01% NP-40, 0.2 mM PMSF, 0.5 mM benzamidine, 1 mM DTT, 15% glycerol, 2 µg/ml leupeptin, 1 µg/ml aprotinin, 0.4 mg/ml recombinant human insulin, 0.4 mg/ml 3x FLAG peptide.

Molarity: 4.7 µM

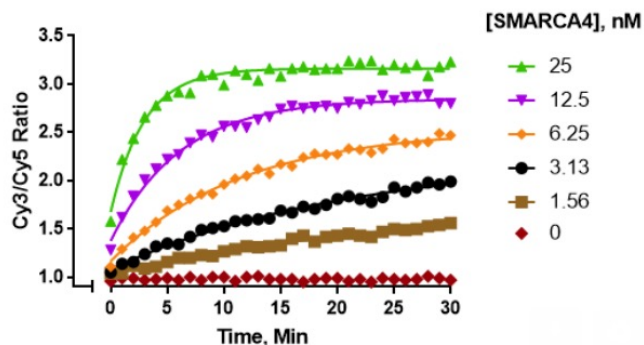
Storage and Stability:

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

Application Notes:

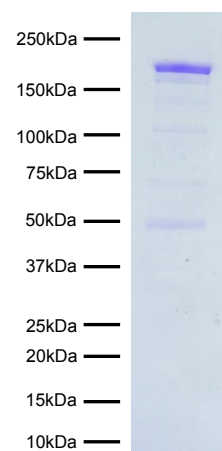
This product is sufficient to perform 100 remodeling reactions using EpiDyne®-FRET substrate (Catalog No. 16-4201). A single reaction is defined as 10 µl containing [20 nM SMARCA4/ 20 nM EpiDyne®-FRET/ 2.5 mM rATP] and remodels to completion in <20 minutes. 5x SMARCA Remodeling Assay Buffer is included (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 best results, make single-use aliquots when first thawed.

References:



ATP-dependent Chromatin Remodeling Assay:

EpiDyne®-FRET Chromatin Remodeling Substrate (20 nM; Catalog No. 16-4201) incubated with SMARCA4 Remodeling Enzyme (concentrations indicated). Curves denote FRET efficiency/chromatin remodeling.



Protein Gel Data: SMARCA4 Remodeling Enzyme (1µg) 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.