

Product Information

Description:	Recombinant human SMARCA2 (SWI/SNF related, matrix associated, actin dependent regulator of chromatin, subfamily a, member 2), encompassing amino acids 636-1331. This construct contains an N-terminal His-tag (6xHis). This protein was affinity purified.
Background:	SMARCA2 (SWI/SNF related, matrix associated, actin dependent regulator of chromatin, subfamily a, member 2) is a member of the helicase family of proteins, involved in regulating the transcription of certain genes and their chromatin structure, DNA repair and genome instability. SMARCA2 genes are often silenced in tumors, resulting in low expression of this protein in most cases. High expression of SMARCA2 has been linked to a good prognosis and to benign differentiated tumors. The selective targeting of SMARCA2 versus SMARCA4 has been a challenge. The use of PROTAC-mediated degradation has proved an alternative for specific degradation of SMARCA2.
Species:	Human
Construct:	SMARCA2 (His-636-1331)
Concentration:	0.35 mg/ml
Expression System:	Sf9
Purity:	83%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	82 kDa
Genbank Accession:	NM_003070
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

