

## Data Sheet

### SMARCA2 (1375-1511)

Human, recombinant, N-terminal HIS-tag  
**Catalog #:** 31111  
Lot #: 120327-GF      Conc.: 0.65 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 0.04 % Tween-  
20, 20% glycerol.

**Stability:** >6 months at -80 °C

#### References:

1. Shain, A.H., *et al.*, *Proc Natl Acad Sci USA*. 2012 Jan 31; **109(5)**: E252-9.
2. Liu, G., *et al.*, *Oncogene*. 2011 Jul 21; **30(29)**: 3295-304.

#### Description:

Human Bromodomain SMARCA2 or BRM, amino-acids 1375 – 1511 (GenBank Accession No. NM\_003070) with N-terminal HIS-tag, MW = 16.9 kDa, expressed in an *E. coli* expression system.

#### Application:

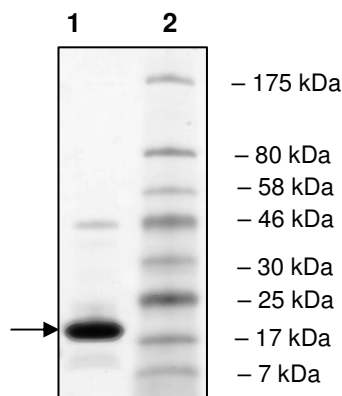
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

## Quality Assurance

### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 µg SMARCA2  
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 16.9 kDa  
**Purity:** ≥92%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MHHHHHKLSPNPPKLTQMNAIIDTVINYKDRCNVEKVP  
SNSQLEIEGNSSGRQLSEVFIQLPSRKELPEYYELIRKPV  
DFKKIKERIRNHKYRSLGDLEKDVMLLCHNAQTFNLEGSQ  
IYEDSIVLQSVFKSARQKIAKEEE

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)