

## Data Sheet

### **SMARCA4/BRG1 (1480-1603)**

Human, recombinant, N-terminal GST tag

**Catalog #:** 31132

Lot #: 121025-D      Conc.: 2.88 mg/ml

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

**Stability:** >6 months at -80°C

**References:**

1. Curtis, C.D., *et al.*, *Mol Cell Biol.* 2012 Apr; **(32)**7: 1312-1320.
2. He H., *et al.*, *Cell Cycle* 2012 Feb 1; **11**(3): 617-627.

**Description:**

Human BRG1, also known as SMARCA4, amino-acids 1480 – 1603 (GenBank Accession No. NM\_1128849) with N-terminal GST tag, MW = 41.3 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:**

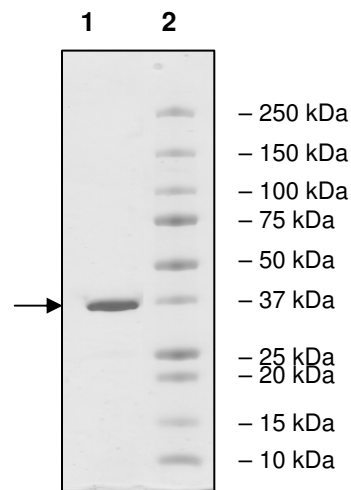
2.5 µg BRG1

**Lane 2:**

Protein Marker

**MW:** 41.3 kDa

**Purity:** ≥96%



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)

**Sequence:**

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMARIYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSDFETLK  
VDFLSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSAEKLS  
PNPPNLTKMKKIVDAVIKYKDSSSGRQLSEVFIQLPSRKELPEYYELIRKPVDFKKIKE  
RIRNHKYRSLNDLEKDVMLLCQNAQTFNLEGLIYEDSIVLQSVFTSVRQKIEKEDDSE

Red: GST tag

Blue: Linker including PreScission Protease Cleavage Site

Black: BRG1 (a.a. 1480-1603)

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)