

## Data Sheet

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

Human, recombinant, N-terminal HIS tag

**Catalog #:** 31102

Lot #: 130411      Conc.: 2.34 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 8.0,  
138 mM NaCl, 2.7 mM KCl, <0.01% Tween-  
20, and 10% 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 HIS tag, MW = 15.5 kDa, expressed in an *E. coli* expression system.

**Application:**

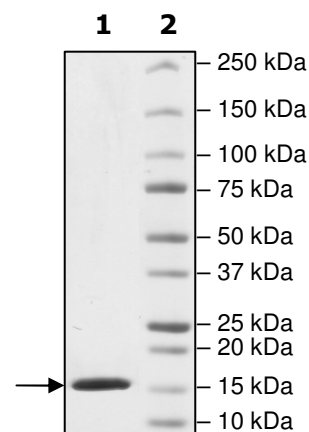
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

## **Quality Assurance**

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

**Lane 1:**  
2 µg BRG1  
**Lane 2:**  
Protein Marker

**MW:** 15.5 kDa  
**Purity:** ≥99%



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)