

## Data Sheet

### SMARCA2, His-Tag, Biotin-Labeled

Human, Recombinant, N-terminal Avi-His-tag

**Catalog #:** 31129

**Lot #:** 160902

**Conc.:** 0.80 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:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended. If necessary, use carrier protein (BSA 0.1-0.5%).

**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:**

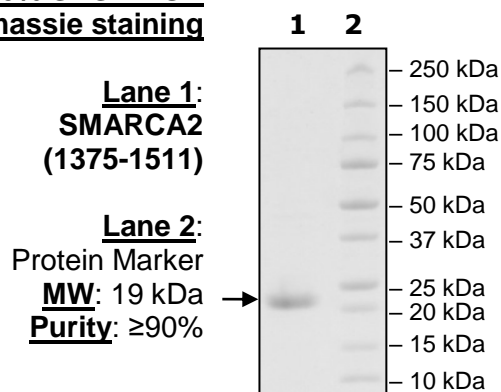
Human Bromodomain SMARCA2 or BRM, a.a. 1375 – 1511 (GenBank Accession No. NM\_003070) with N-terminal Avi-His-tag, MW = 19 kDa, expressed in an *E. coli* expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ .

**Application:**

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

## **Quality Assurance**

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



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)