

Data Sheet

BRPF1 (627-746)

Human, recombinant, N-terminal HIS-tag

Catalog #: 31112

Lot #: 130412 Conc.: 2.18 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, <0.01 % Tween-
20, and 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Laue, K., *et al.*, *Development*. 2008
June; **(135)111**: 1935-46.
2. Mayya, V., *et al.*, *Sci Signal*. 2009 Aug
18; **2(84)**: ra46.

Description:

Human Bromodomain and PHD finger
containing 1, or BRPF1, amino-acids 627
– 746 (GenBank Accession No.
NM_001003694) with N-terminal HIS-tag,
MW = 15.1 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:

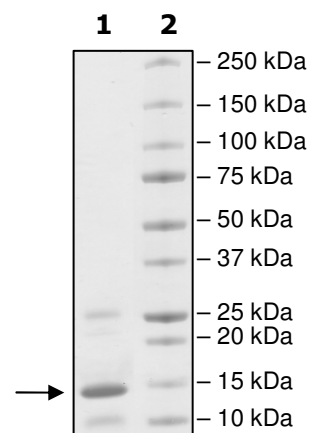
2 µg BRPF1

Lane 2:

Protein Marker

MW: 15 kDa

Purity: ≥68%



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

Please visit our website at: 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

Protein Sequence:

MHHHHHMQLTPFLILLRKTLEQLQEKDTGNIFSEPVPLSEVTELDEVPDYLDHIKKPMD
FFTMKQNLEAYRYLNFDDFEEDFNLIVSNCLKYNAKDTIFYRAAVRLREQGGAVLRQAR
RQAEKMG

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
Please visit our website at: www.bpsbioscience.com