

Data Sheet

BRPF3 (576-701)

Human, recombinant, N-terminal GST-tag

Catalog #: 31130

Lot #: 130326-G1 Conc.: 0.39 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. Ullah, M., *et al.*, *Mol Cell Biol.* 2008 Nov;**28(22)**:6828-6843.
2. Mishima, Y., *et al.*, *Blood.* 2011 Sep 1;**118(9)**:2443-53.
3. Slama, P., *et al.*, *Nucleic Acids Res.* 2011 Mar;**39(5)**:1666-1679.

Description:

Human Bromodomain and PHD finger containing 3, or BRPF3, amino-acids 576 – 701 (GenBank Accession No. NM_015695) with N-terminal GST-tag, MW = 41.2 kDa, expressed in an *E. coli* expression system.

Application:

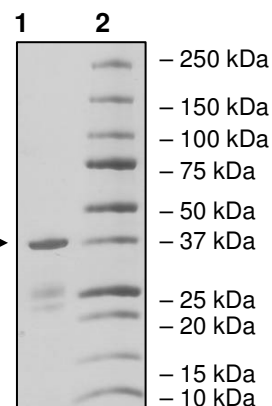
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg BRPF3
Lane 2:
Protein Marker

MW: 41.2 kDa →
Purity: ≥80%



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

Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQATFGGGDHPPKSDLEVLFQGPLGSQVKV
QQAAMELELMPFNVLLRRTLDLLQEKP AHIFAEPVNLSEVPDYLEFISKPMDFSTMR
RKLESHLYRTLEEFEEFNLIVTNCMKYNAKDTIFHRAAVRLRDLGGAILRHARRQAE
NIG

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