

Data Sheet

BRD3, BD1 and BD2 (29-417)

Human, recombinant, N-terminal GST tag

Catalog #: 31035

Lot #: 120810-G2D Conc.: 0.18 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:

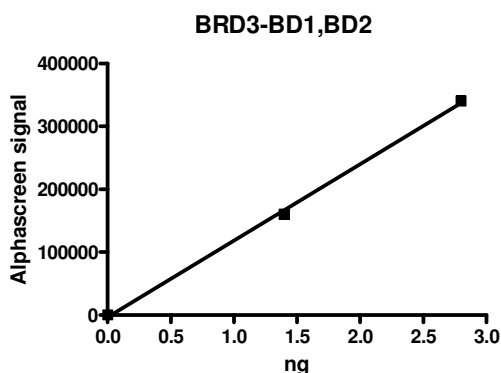
1. Gamsjaeger, R., *et al.*, *Mol Cell Biol.* 2011 Jul; **31(13)**: 2632-2640.
2. Lamonica, JM, *et al.*, *Proc Natl Acad Sci USA.* 2011 May 31; **108(22)**: E159-168.

Description: Human Bromodomain containing 3, also known as BRD3 or RING3L, amino-acids 29 – 417 (GenBank Accession No. NM_007371) with N-terminal GST tag, MW = 70 kDa, expressed in an *E. coli* expression system.

Assay Conditions: BRD3 protein is incubated with biotinylated acetyl-histone peptide substrate (Cat. #33000) in BRD assay buffer (Cat. #33001) in a 10 µl reaction for 30 minutes at RT. AlphaLISA® GSH acceptor beads (Perkin Elmer) are added, followed by AlphaScreen® Streptavidin-conjugated donor beads (Perkin Elmer). After 1 hour incubation at RT, Alphacounts are measured.

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

Quality Assurance

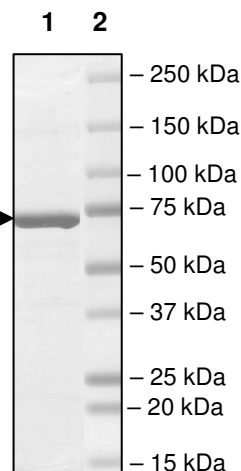


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 µg BRD3

Lane 2:
 Protein Marker
 Biorad (#161-0373)

MW: 70 kDa
Purity: ≥77%



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

Amino Acid Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPOIDKYLKSSKYIAWPLQGQATFGGGDHPPKSDLEVLFGGPLGSPSKPG
RKTNQLQYMQNVVVKTLWKHQFAWPFYQPVDAIKLNLDPYHKIKNPMDMGTIKKRL
ENNYYSASECMQDFNTMFTNCYIYNKPTDDIVLMAQALEKIFLQKVAQMPQEEVEL
LPPAPKGKGRKPAAGAQSAGTQQVAAVSSVSPATPFQSVPTVSQTPVIAATPVPTIT
ANVTSVPVPPAAAPPPATPIVPVVPPTPPVVKKKGVKRKADTTTPTTSAITASRSESP
PPLSDPKQAKVVARRESGGRPIKPPKDLLEDGEVPQHAGKKGKLEHLRYCDSILRE
MLSKKHAAYAWPFYKPVDAEALHDYHDIKHMPMDLSTVKKRKMMDGREYPDAQGFA
ADVRLMFSNCYKYNPPDHEVVAMARKLQDVFEMRFAKMPD

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