

Data Sheet

BRD4, BD1 and BD2 (49-460)

Human, recombinant, N-terminal His tag

Catalog #: 31045

Lot #: 120524-G Conc.: 0.34 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8,
 124 mM NaCl, 2.4 mM KCl, and 10%
 glycerol.

Stability: >6 months at -80 °C

References:

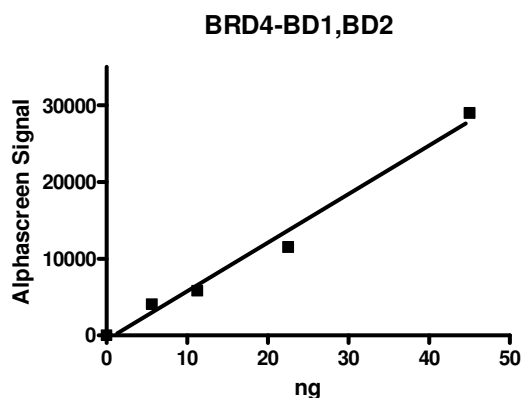
- Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012 May; **90(5)**: 587-95.
- Voigt, P, *et al.*, *Genome Biol*. 2011 Nov 29; **12(11)**:133.

Description: Human Bromodomain containing 4, also known as BRD4, HUNK1, or MCAP, amino-acids 49-460 (GenBank Accession No. NM_058243) with N-terminal His tag, MW = 47.5 kDa, expressed in an *E. coli* expression system.

Assay Conditions: BRD4 protein is incubated with biotinylated acetyl-histone peptide substrate (Cat. #33000) in BRD assay buffer (Cat. #33001) in a 10 µl reaction for 1 hour at RT. AlphaLISA® Nickel Chelate 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.

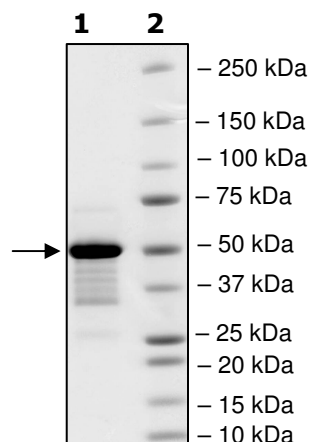
Quality Assurance



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

Lane 1:
 4 µg BRD4
Lane 2:
 Protein Marker
 Biorad (#161-0373)

MW: 47.5 kDa
Purity: ≥75%



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

MHHHHHHETSNPNKPKRQTNQLQYLLRVVLKTLWKHQFAWPFQQPVDAVKLNLPD
YYKIIKTPMDMGTIKKRENNYYWNAQECIQDFNTMFTNCYIYNKPGDDIVLMAEALEK
LFLQKINELPTEETEIMIVQAKGRGRGRKGTAKPGVSTVPNTTQASTPPQTQTPQP
NPPPQATPHPPFAVTPDLIVQTPVMTVVPPQPLQTPPPVPPQPQPPPAPAPQPVS
HPPIAATPQPVKTKKGVKRKADTTTPTTIDPIHEPPSLPPEPKTTKLGQRRESSRPVK
PPKDVPSQQHPAPEKSSKVSEQLKCCSGILKEMFAKKHAAYAWPFYKPVDEAL
GLHDYCDIHKPMDMSTIKSKLEAREYRDAQEFGADVRLMFSNCYKYNPPDHEVVAM
ARKLQDVFEMRFAKMPDE

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