

Data Sheet

BRD4, BD1 and BD2 (49-460)

Human, recombinant, un-tagged
 Catalog #: 31047
 Lot #: 130815-M Conc.: 3.0 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol.

Stability: >6 months at -70°C. Avoid multiple freeze/thaw cycles.

References:

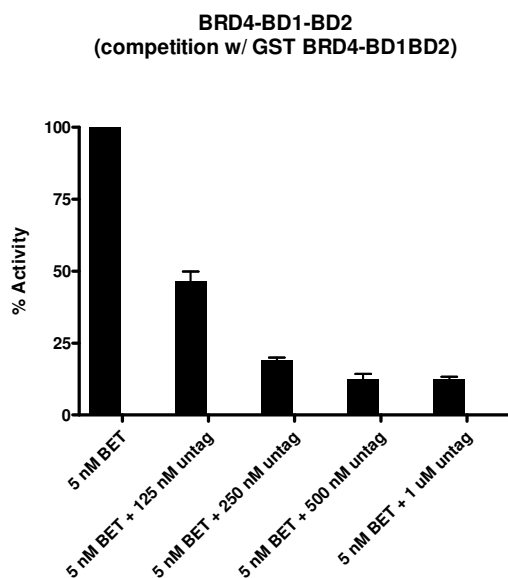
- Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012 May; **90(5)**: 587-95.
- Voigt, P, *et al.*, *Genome Biol*. 2011 Nov 29;**12(11)**:133.

Description: Human Bromodomain containing 4, also known as BRD4, HUNK1, or MCAP, amino-acids 49-460 (GenBank Accession No. NM_058243) with no tag, MW = 46.9 kDa, expressed in an *E. coli* expression system.

Assay Conditions: BET Ligand (Cat. # 33000) was incubated with un-tagged BRD4 for 15 min at room temperature in 1X BRD assay buffer. GST-tagged BRD4 (Cat.# 31044) was added and incubated for 20 min followed by addition of 20 µl mixture of GSH-Acceptor beads and Streptavidin-Donor beads for AlphaScreen® analysis. Total reaction mixture was incubated for 35 min and the AlphaScreen® signal was measured by EnSpire® Multilabel Plate Reader (PerkinElmer).

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

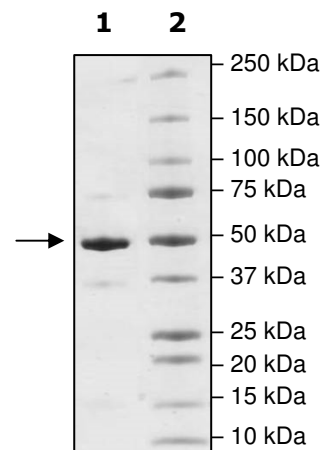
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg BRD4
Lane 2:
Protein Marker

MW: 46.9 kDa
Purity: ≥90%



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

**GPLGSETSNPNKPKRQTNQLQYLLRVVLKTLWKHQFAWPFQQPVDAVKLNLPDYK
IIKTPMDMGTIKKRENNYYWNAQECIQDFNTMFTNCYIYNKPGDDIVLMAEALEKFL
QKINELPTEETEIMIVQAKGRGRGRKETGTAKPGVSTVPNTTQASTPPQTQTPQPNPP
PVQATPHFPFAVTPDLIVQTPVMTVVPPQPLQTPPPVPPQPQPPAPAPQPVSHPPII
AATPQPVKTKKGVKRKADTTTPTTIDPIHEPPSLPPEPKTTKLGQRRESSRPVKPPK
DVPDSQQHPAPEKSSKVSEQLKCCSGILKEMFAKKHAAYAWPFYKPDVEALGLHD
YCDIIKHPMDMSTIKSKLEAREYRDAQEFGADVRLMFSNCYKYNPPDHEVVAMARKL
QDVFEMRFAKMPDE**

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