

Data Sheet

BRD4 (342-460)

Human, recombinant, N-terminal GST tag

Catalog #: 31041

Lot #: 121109-D Conc.: 3.7 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:

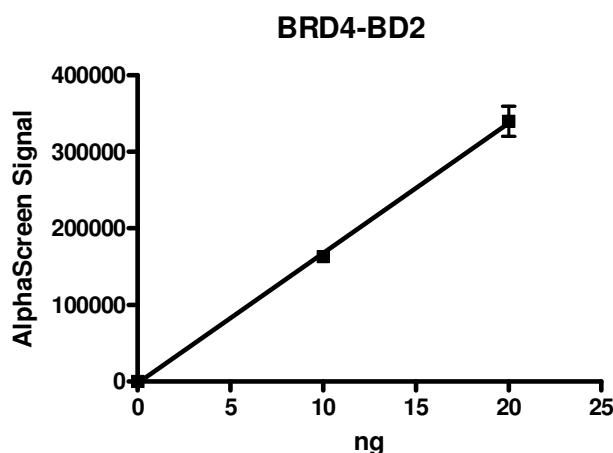
1. Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012 May; **90(5)**: 587-95.
2. Voigt, P, *et al.*, *Genome Biol.* 2011 Nov 29; **12(11)**:133.

Description: Human Bromodomain containing 4, also known as BRD4, HUNK1, or MCAP, amino-acids 342-460 (GenBank Accession No. NM_058243) with N-terminal GST tag, MW = 40.6 kDa, expressed in an *E. coli* expression system.

Assay Conditions: Assay was performed using BRD4 Inhibitor Screening Assay Kit, Cat. #32524. 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® GSH acceptor beads (Perkin Elmer) are added, followed by AlphaScreen® Streptavidin-conjugated donor beads (Perkin Elmer). After 15 minute incubation at RT, Alphacounts are measured.

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

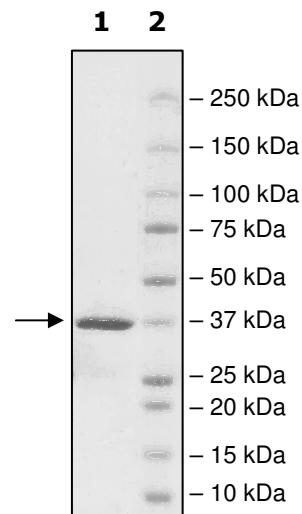
Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:
4 µg BRD4
Lane 2:
Protein Marker

MW: 40.6 kDa
Purity: ≥96%



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