

Data Sheet

BRD4, BD1 and BD2 (49-460)

Human, recombinant, N-terminal GST tag

Catalog #: 31044

Lot #: 140529-2

Conc.: 0.34 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 -80°C. *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

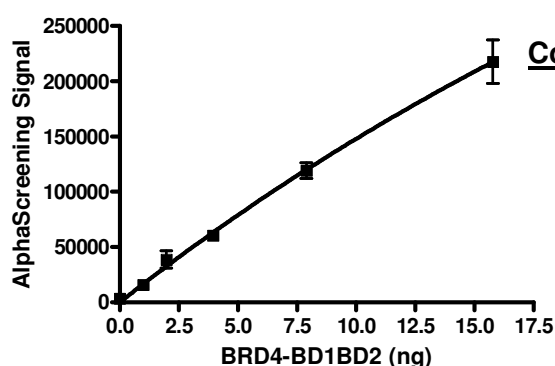
1. Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012 May;**90(5)**:587-95.
2. Voigt, P, *et al.*, *Genome Biol*. 2011 Nov 29;**12(11)**:133.

Description: Human Bromodomain containing 4, also known as BRD4, HUNK1, or MCAP, amino-acids 49-460 (GenBank Accession No. NM_058243) with N-terminal GST tag, MW = 73.3 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® 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.

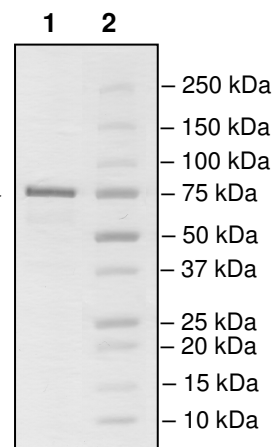
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 3 µg BRD4
Lane 2: →
 Protein Marker

MW: 73.3 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