

## Data Sheet

### BRD4 (BD2), His-tag

Human, recombinant, N-terminal His-tag

**Catalog #:** 31043

**Lot #:** 190509

**Conc.:** 2.38 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 8.0, 138 mM NaCl, 2.7 mM KCl, 0.05% Tween-20, and 10% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

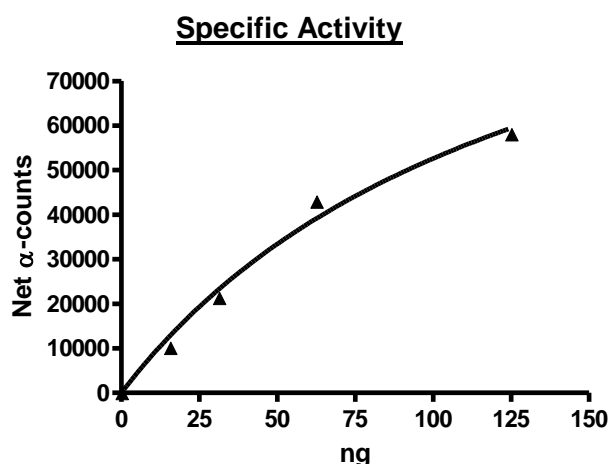
- Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012 May; **90(5)**:587-95.
- Voigt, P, *et al.*, *Genome Biol*. 2011 Nov 29; **12(11)**:133.

**Description:** Human BRD4 also known as Human Bromodomain containing 4, HUNK1, and MCAP, GenBank Accession No. NM\_058243, a.a. 342-460 with N-terminal His-tag, expressed in an *E. coli* expression system. MW = 15 kDa.

**Assay Conditions:** BRD4 protein is incubated with biotinylated acetyl-histone peptide substrate 2.5  $\mu\text{M}$  stock (BPS Bioscience #33000) in BRD assay buffer (BPS Bioscience #33001) in a 10  $\mu\text{l}$  reaction for 1 hour at room temperature. AlphaLISA<sup>®</sup> Nickle acceptor beads (PerkinElmer) are added, followed by Alphascreen<sup>®</sup> Streptavidin-conjugated donor beads (PerkinElmer). After 1 hour of incubation at room temperature, Alphacounts are measured.

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

## Quality Assurance

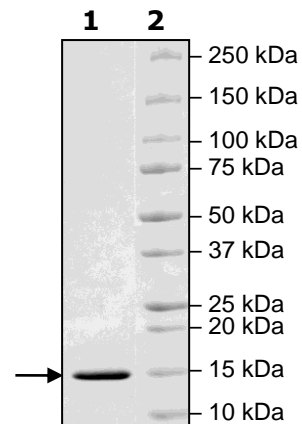


**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
2  $\mu\text{g}$  BDR4(BD2)

**Lane 2:**  
Protein Marker

**MW:** 15 kDa  
**Purity:**  $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)