

## Data Sheet

### **BRD4 (49-170)**

Human, recombinant, N-terminal His-tag

**Catalog #:** 31042

**Lot #:** 140918-2      **Conc.:** 5.39 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  $-80^{\circ}\text{C}$ . *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein BSA (0.1 – 0.5%).*

#### References:

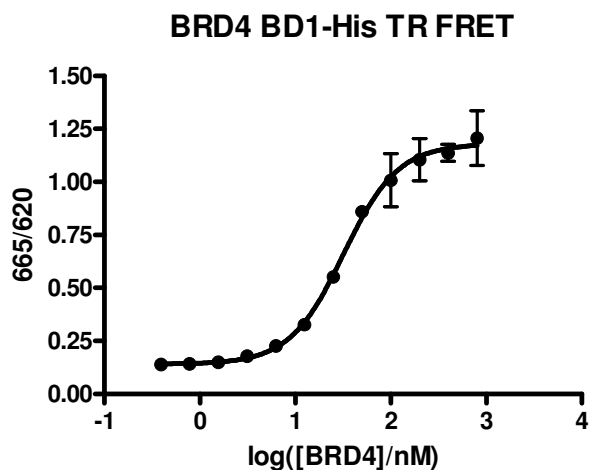
1. Rodriguez, RM, *et al.*, *J Mol Med (Berl)*. 2012; **90(5)**: 587-95.
2. Voigt, P, *et al.*, *Genome Biol*. 2011; **12(11)**:133.

**Description:** Human Bromodomain containing 4, also known as BRD4, HUNK1, or MCAP, amino-acids 49-170 (GenBank Accession No. NM\_058243) with N-terminal His tag, MW = 15.5 kDa, expressed in an *E. coli* expression system.

**Assay Conditions:** Varying concentrations of BRD4 are incubated with 200 nM BET Bromodomain Ligand, Tb donor and dye-labeled acceptor, in BRD TR-FRET assay buffer, in a 20  $\mu\text{l}$  reaction for 2 hours at RT. Donor emission at 620 nm followed by dye-acceptor emission at 665 nm are measured and analysis is performed using the 665/620 ratio.

**Application:** Useful for the study of bromodomain binding assays, screening

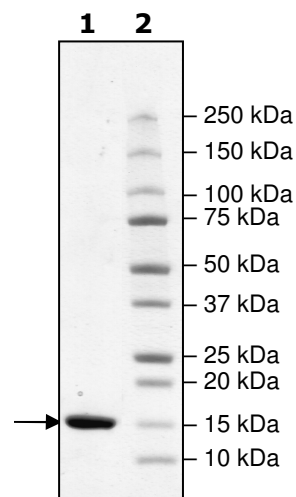
## Quality Assurance



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

**Lane 1:**  
2  $\mu\text{g}$  BRD4  
**Lane 2:**  
Protein Marker

**MW:** 15.5 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 **1.858.829.3082** 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)