

## Data Sheet

### **BRD3 (29-145)**

Human, recombinant, N-terminal HIS tag

**Catalog #:** 31030

**Lot #:** 120207G      **Conc.:** 0.52 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:** >6 months at -80 °C

#### **References:**

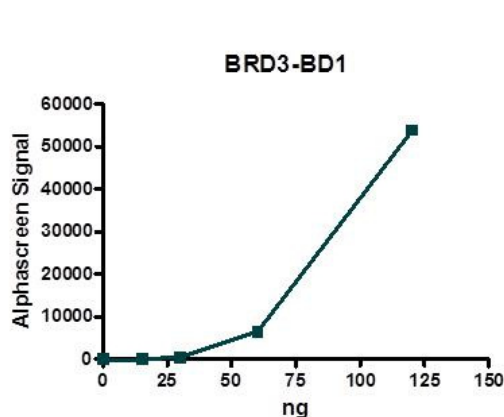
1. Gamsjaeger, R., *et al.*, *Mol Cell Biol.* 2011 Jul; **31(13)**: 2632-2640.
2. Lamonica, JM, *et al.*, *Proc Natl Acad Sci USA.* 2011 May 31; **108(22)**: E159-168.

**Description:** Human Bromodomain containing 3, also known as BRD3 or RING3L, amino-acids 29-145 (GenBank Accession No. NM\_007371) with N-terminal HIS tag, MW = 14.9 kDa, expressed in an *E. coli* expression system.

**Assay Conditions:** BRD3 is incubated with biotinylated acetyl-histone peptide substrate (BPS #33000) in BRD assay buffer (BPS #33001) in a 10 µl reaction for 1 hour at RT. AlphaLISA® Nickel Chelate 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**



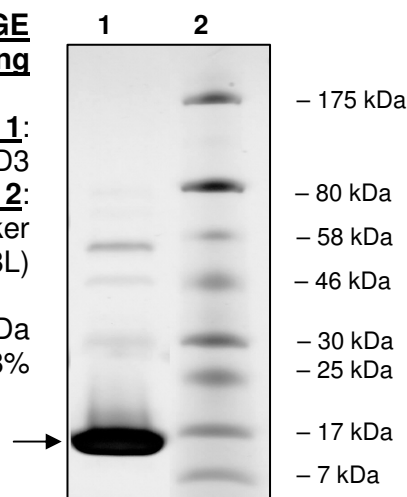
#### **4-20% SDS-PAGE** **Coomassie staining**

**Lane 1:**  
4.5 µg BRD3

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 14.9 kDa

**Purity:** ≥88%



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)