

## Data Sheet

### BRD9 (135-242)

Human, recombinant, N-terminal GST tag  
**Catalog #:** 31091  
**Lot#:** 130315-D      **Conc.:** 3.6 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. Avoid multiple freeze/thaw cycles.

**References:**

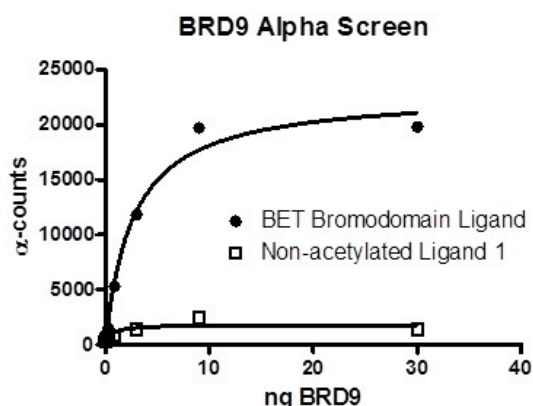
1. Scotto, L., *et al.*, *Mol. Cancer*. 2008 Jun 17; **7**:58.
2. Kang, J.U., *et al.*, *Cancer Genet Cytogenet*. 2008 Apr 1; **182(1)**: 1-11.

**Description:** Human Bromodomain containing 9, also known as BRD9, amino-acids 135 – 242 (GenBank Accession No. NM\_023924) with N-terminal GST tag, MW = 39 kDa, expressed in an *E. coli* expression system.

**Assay Conditions:** Assay was performed using BRD9 Inhibitor Screening Assay Kit, Cat. # 32519. BRD9 protein is incubated with biotinylated acetyl-histone peptide substrate BET Bromodomain Ligand in BRD9 assay buffer in a 10 ul reaction for 30 min at RT. AlphaLISA® GSH acceptor beads (PerkinElmer) were then added, and the mixture was incubated for 30 minutes at room temperature (RT). AlphaScreen® Streptavidin-conjugated donor beads (PerkinElmer) were then added. After 10 minute incubation at RT, Alphacounts are measured.

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

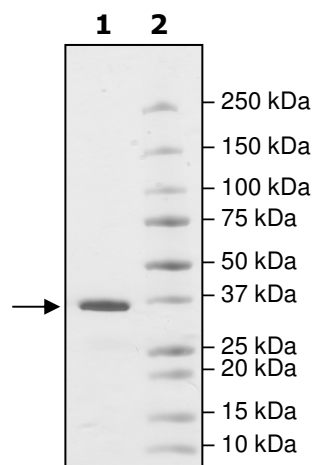
## Quality Assurance



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

**Lane 1:**  
 2 µg BRD9  
**Lane 2:**  
 Protein Marker

**MW:** 39 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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYID  
GDVKLTQSMARIIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLKVDF  
LSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLVCFKK  
RIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDLEVLFGGPLGSENESTPIQQLL  
EHFLRQLQRKDPHGFFAFPVTDIAIPGYSMIIKHPMDFGTMKDKIVANEYKSVTEFKADF  
KLMCDNAMTYNRPDVTYYYKLAKKILHAGFKMMSKQAA

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)