

Data Sheet

BRD9 (135-242)

Human, recombinant, N-terminal HIS tag

Catalog #: 31090

Lot #: 140218

Conc.: 1.7 mg/ml

Formulated in: 40 mM Tris-HCl, 110 mM NaCl, 2.2 mM KCl, 0.02% Tween-20 and 20% glycerol.**Stability:** >6 months at -80°C. *Avoid freeze/thaw cycles.***References:**

1. Scotto, L., *et al.*, *Mol. Cancer*. 2008 Jun 17; **7**:58.
2. Kang, JU., *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 HIS tag, MW = 13.4 kDa, expressed in an *E. coli* expression system.**Application:**

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

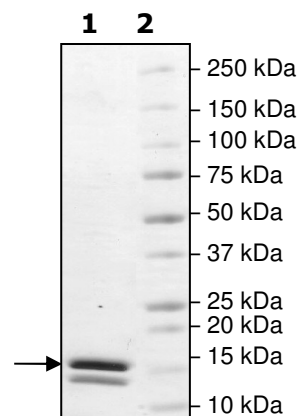
Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:**

2 µg BRD9

Lane 2:

Protein Marker

MW: 13.4 kDa**Purity:** ≥72%

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.comPlease visit our website at: www.bpsbioscience.com