

## Data Sheet

### BRDT, BD1 and BD2 (22-382)

Human, recombinant, N-terminal GST tag  
**Catalog #:** 31061  
Lot #: 121012-GC      Conc.: 0.36 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

**References:**

1. Pivot-Pajot C., *et al.*, *Mol Cell Biol.* 2003 Aug; **23(15)**: 5354-65.
2. Morinière J., *et al.*, *Nature.* 2009 Oct 1; **461(7264)**:664-8.

**Description:**

Human Bromodomain testis-specific protein, also known as BRDT, BRD6, and CT9, amino-acids 22 – 382 (GenBank Accession No. NM\_207189) with N-terminal GST tag, MW = 68.1 kDa, expressed in an *E. coli* expression system.

**Application:**

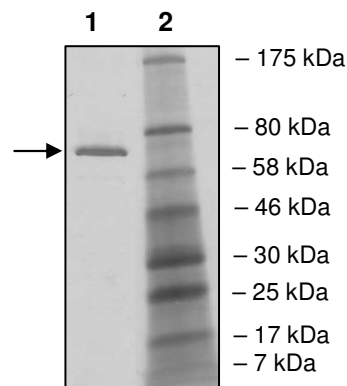
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

## Quality Assurance

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

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

**MW:** 68.1 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:**

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK  
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSTKKN  
RLTNQLQYLQKVVLKDLWKHSFSWPFQRPVDAVKLQLPDYYTIIKNPMDLNTIKRLE  
NKYYAKASECIEDFNTMFSNCYLYNKPGDDIVLMAQALEKLFMQKLSQMPQEEQVVG  
VKERIKKGTQQNIAVSSAKEKSSPSATEKVFKQQEIPSVFPKTSISPLNVVQGASVNSS  
SQTAAQVTKGVKRAKADTTTPATSAVKASSEFSPTFTEKSVLPPIKENMPKNVLPDS  
QQQYNVVKTVKVTEQLRHCSEILKEMLAKKHFSYAWPFYNPVDVNALGLHNYVDVVK  
NPMDLGTIKEKMDNQEYKDAYKFAADVRLMFMNCYKYNPPDHEVVTMARMLQDVFE  
THFSKIPIEPVE

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)