

Data Sheet

BRDT (257-382)

Human, recombinant, N-terminal GST tag

Catalog #: 31060

Lot #: 121011-G Conc.: 0.32 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 257 – 382 (GenBank Accession No. NM_207189) with N-terminal GST tag, MW = 41.7 kDa, expressed in an *E. coli* expression system.

Application:

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

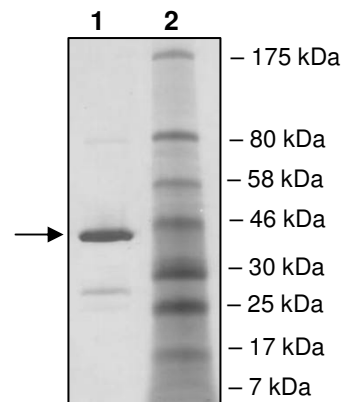
Quality Assurance

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

Lane 1:
2 µg BRDT

Lane 2:
Protein Marker

MW: 41.7 kDa
Purity: ≥85%



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

Please visit our website at: 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

Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSSQQQ
YNNVKTVKVTEQLRHCSEILKEMLAKKHFSYAWPFYNPVDVNALGLHNYDVVKNPM
DLGTIKEKMDNQEYKDAYKFAADVRLMFMNCYKYNPPDHEVVTMARMMLQDVFETHF
SKIPIEPVE

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
Please visit our website at: www.bpsbioscience.com