

## Data Sheet

### BRDT (22-138)

Human, recombinant, N-terminal GST tag

**Catalog #:** 31108

Lot #: 130117-D      Conc.: 3.1 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
 110 mM NaCl, 2.2 mM KCl, 20% Glycerol.

**Stability:** >6 months at -80°C

**References:**

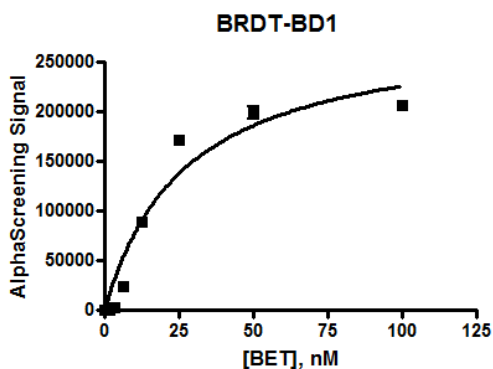
1. Pivot-Pajot C., *et al.*, *Mol Cell Biol.* 2003 Aug; **23(15)**: 5354-65.
2. Morinière J., *et al.*, *Nature.* 2009 Oct 1; **461(7264)**:664-8.

**Description:** Human Bromodomain testis-specific protein, also known as BRDT, BRD6, and CT9, amino-acids 22 – 138 (GenBank Accession No. NM\_207189) with N-terminal GST tag, MW = 40.7 kDa, expressed in an *E. coli* expression system.

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

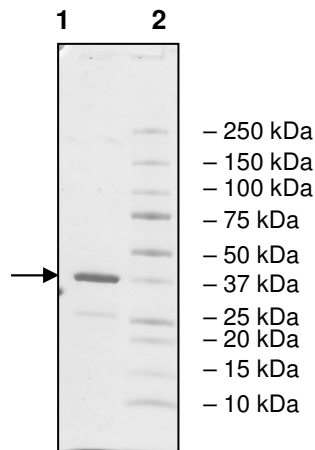
## Quality Assurance



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

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

**MW:** 40.7 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)