

Data Sheet

BRDT (257-382)

Human, recombinant, N-terminal HIS tag

Catalog #: 31100

Lot #: 120216G Conc.: 0.21 mg/ml

Formulated in: 50 mM Tris-HCl, pH 8.0,
138 mM NaCl, 2.7 mM KCl, 0.05% Tween-
20, and 10% 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 HIS tag, MW = 15.8 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:

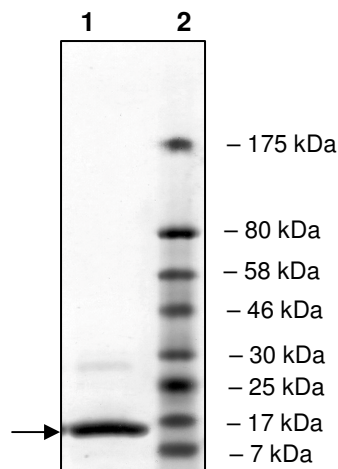
4.2 µg BRDT

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 15.8 kDa

Purity: ≥97%



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