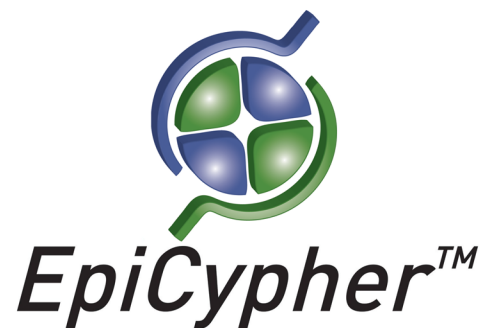


## BRDT Bromodomain 2

**Catalog No.** 15-0017  
**Lot No.** 13228001  
**Pack Size** 100 µg



**Type** Bromodom **Expressed In** *E. coli*  
**Mol. Wgt.** 38 kDa **Epitope Tag** GST

### Product Description:

Recombinant human BRDT bromodomain 2 (CT9, Fsg3, accession number Q58F21, amino acids 258 to 387) containing an N-terminal GST tag, expressed in *E. coli*. BRDT is a testis-specific bromodomain protein that binds histone H4 acetylated at lysine 5 and lysine 8. BRDT is required for spermatogenesis and male mice lacking BRDT are sterile. BRDT is expressed in a percentage of cancers such as non-small cell lung cancer and squamous cell carcinomas of the head, neck and esophagus.

### Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

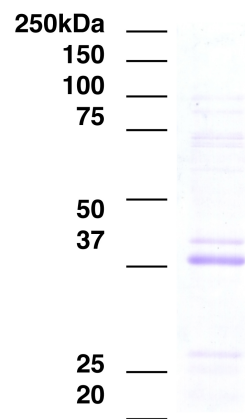
### Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

### Application Notes:

Recombinant BRDT Bromodomain 2 is useful for protein binding and screening experiments examining acetylated protein substrates.

### References:



**Protein Gel Data:** BRDT Bromodomain 2 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.