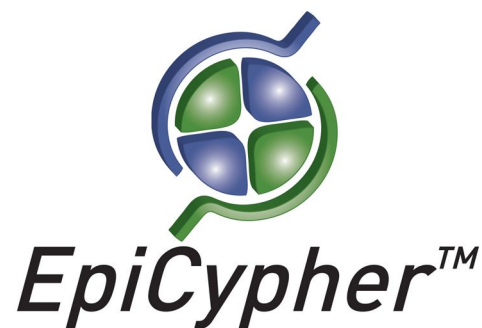


# BRD7 Bromodomain, Recombinant Human

**Catalog No.** 15-0067  
**Lot No.** 17041001  
**Pack Size** 100 µg



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

## Product Description:

BRD7 Bromodomain, Recombinant Human (Bromodomain- containing protein 7, BP75, CELTIX-1, accession number Q9NPI1, amino acids 127 - 241) containing an N-terminal GST tag, expressed in *E. coli*. BRD7 is part of the PBAF SWI/SNF chromatin remodeling complex. BRD7 bromodomain interacts with acetylated histone H3 (lys 9 and lys 14) and H4 (lys 8). BRD7 has been reported to function as both a co-activator and a co-repressor of transcription.

## Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 50 mM TRIS pH 7.5, 200 mM NaCl, 1 mM DTT, 20% glycerol.

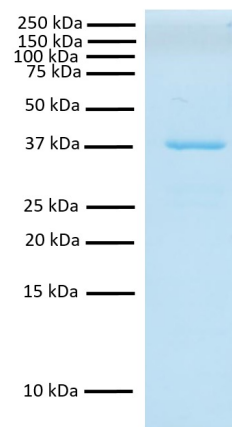
## Storage and Stability:

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

## Application Notes:

BRD7 Bromodomain, Recombinant Human is useful for protein binding and screening experiments examining acetylated protein and peptide substrates.

## References:



**Protein Gel Data:** BRD7 Bromodomain, Recombinant Human (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.