

# PRODUCT INFORMATION



## BRD2 bromodomain 2 (human, recombinant)

Item No. 11070

### Overview and Properties

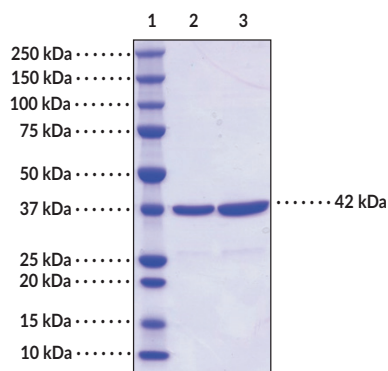
**Synonyms:** Bromodomain containing protein 2, RING3, RNF3  
**Source:** Recombinant N-terminal GST-tagged protein expressed in *E. coli*  
**Amino Acids:** 339-459  
**Uniprot No.:** P25440  
**Molecular Weight:** 42 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥2 years  
**Purity:** *batch specific* (≥95% estimated by SDS-PAGE)  
**Supplied in:** 50 mM Tris, pH 7.5, with 500 mM sodium chloride, 5% glycerol, and 5 mM β-mercaptoethanol

### Protein

**Concentration:** *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: BRD2 Domain 2 (2 μg)  
Lane 3: BRD2 Domain 2 (4 μg)

Representative gel image shown; actual purity may vary between each batch.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/16/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

Bromodomain-containing 2 (BRD2) is a transcriptional regulator that is a member of the bromodomain and extra-terminal (BET) family.<sup>1</sup> It is ubiquitously expressed and localizes to the nucleus. BRD2 is composed of two N-terminal bromodomains (BD1 and BD2) that bind acetylated lysine on histones, serving to couple histone acetylation marks to the transcriptional regulation of target promoters, and an extra-terminal domain that mediates chromatin interactions.<sup>1,2</sup> BRD2 associates with transcription effector and regulator proteins, including RNA polymerase II, histone acetylases and deacetylases, and transcriptional co-activators and co-repressors to form a transcription complex that regulates the expression of genes involved in inflammation and cell proliferation.<sup>1,3</sup> BRD2 also binds to the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) envelope (E) protein, a transmembrane protein involved in CoV virion assembly and pathogenesis of the related virus, SARS-CoV.<sup>4-6</sup> The BD2 bromodomain of BRD2 contains four  $\alpha$ -helices that form the binding site for the acetylated histone H4 tail.<sup>7</sup> Point mutations of valine 329 (V329A) or asparagine (N382A) in the BD2 bromodomain eliminate BRD2 binding to acetylated histone H4 lysine 12 (H4K12Ac), and a point mutation of tyrosine 427 (Y427F) abolishes the association between BRD2 and STAT3.<sup>8,9</sup> Cayman's BRD2 bromodomain 2 (human, recombinant) protein can be used for ELISA, Western blot (WB), and binding assay applications.

## References

1. Taniguchi, Y. The bromodomain and extra-terminal domain (BET) family: Functional anatomy of BET paralogous proteins. *Int. J. Mol. Sci.* **17**(11), E1849 (2016).
2. Hnilicová, J., Hozeifi, S., Stejskalová, E., *et al.* The C-terminal domain of Brd2 is important for chromatin interaction and regulation of transcription and alternative splicing. *Mol. Biol. Cell* **24**(22), 3557-3568 (2013).
3. Gilan, O., Rioja, I., Knezevic, K., *et al.* Selective targeting of BD1 and BD2 of the BET proteins in cancer and immuno-inflammation. *Science* (2020).
4. Gordon, D.E., Jang, G.M., Bouhaddou, M., *et al.* A SARS-CoV-2-human protein-protein interaction map reveals drug targets and potential drug-repurposing. *BioRxiv* (2020).
5. Kandeel, M., Ibrahim, A., Fayez, M., *et al.* From SARS and MERS CoVs to SARS-CoV-2: Moving toward more biased codon usage in viral structural and nonstructural genes. *J. Med. Virol.* (2020).
6. Schoeman, D. and Fielding, B.C. Coronavirus envelope protein: Current knowledge. *Viol. J.* **16**(1), 69 (2019).
7. Umehara, T., Nakamura, Y., Wakamori, M., *et al.* Structural implications for K5/K12-di-acetylated histone H4 recognition by the second bromodomain of BRD2. *FEBS Letters* **584**(18), 3901-3908 (2010).
8. Huang, H., Zhang, J., Shen, W., *et al.* Solution structure of the second bromodomain of Brd2 and its specific interaction with acetylated histone tails. *BMC Struct. Biol.* **7**, 57 (2007).
9. Cheung, K.L., Zhang, F., Jaganathan, A., *et al.* Distinct roles of Brd2 and Brd4 in potentiating the transcriptional program for Th17 cell differentiation. *Mol. Cell* **65**(6), 1068-1080 (2017).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM