

PRODUCT INFORMATION



BRD2 bromodomain 1 and 2 (human, recombinant)

Item No. 11069

Overview and Properties

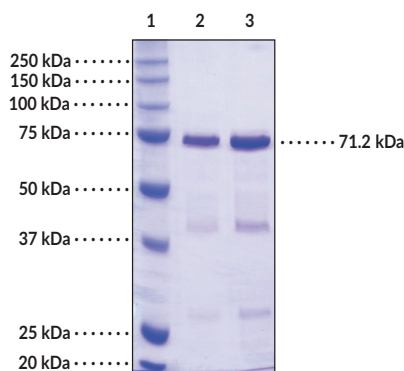
Synonyms: Bromodomain containing protein 2, RING3, RNF3
Source: Recombinant N-terminal GST-tagged protein expressed in *E. coli*
Amino Acids: 65-459
Uniprot No.: P25440
Molecular Weight: 71.2 kDa
Storage: -80°C (as supplied)
Stability: ≥2 years
Purity: *batch specific* (≥65% estimated by SDS-PAGE)
Supplied 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 Domains 1 and 2 (2 μg)

Lane 3: BRD2 Domains 1 and 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, 08/06/2020

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.¹ 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.^{1,2} 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.^{1,3} 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.⁴⁻⁶ siRNA knockdown of *Brd2* reduces LPS-induced production of TNF and IL-6 in mouse bone marrow-derived macrophages (BMDMs).⁷ Genetic deletion of *Brd2* in mice induces neural tube defects and is embryonic lethal.¹ *Brd2*^{+/-} mice have decreased numbers of GABAergic neurons and a reduced flurothyl-induced seizure threshold. SNPs in *BRD2* have been found in individuals with juvenile myoclonic epilepsy. Cayman's BRD2 bromodomains 1 and 2 (human, recombinant) protein can be used for Western blot (WB) 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. *Virol. J.* **16(1)**, 69 (2019).
7. Belkina, A.C., Nikolajczyk, B.S., and Denis, G.V. BET protein function is required for inflammation: Brd2 genetic disruption and BET inhibitor JQ1 impair mouse macrophage inflammatory responses. *J. Immunol.* **190(7)**, 3670-3678 (2013).

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