

# PRODUCT INFORMATION



## BRD4 bromodomains 1 and 2 (human, recombinant; aa 2-477)

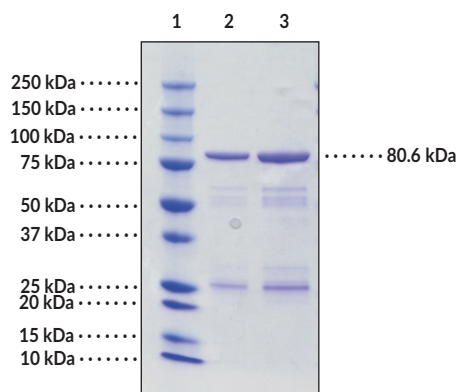
Item No. 14822

### Overview and Properties

<b>Synonyms:</b>	Bromodomain containing protein 4, HUNK1, MCAP, Mitotic Chromosome-associated Protein
<b>Source:</b>	Recombinant N-terminal GST-tagged protein expressed in <i>E. Coli</i>
<b>Amino Acids:</b>	2-477 (partial protein)
<b>Uniprot No.:</b>	O60885
<b>Molecular Weight:</b>	80.6 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥2 years
<b>Purity:</b>	<i>batch specific</i> (≥70% estimated by SDS-PAGE)
<b>Supplied in:</b>	10 mM HEPES, pH 7.5, with 500 mM sodium chloride, and 5% glycerol
<b>Protein Concentration:</b>	<i>batch specific</i> 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: BRD4 Domains 1 and 2 (2 µg)

Lane 3: BRD4 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, 04/29/2020

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 protein 4 (BRD4) is a member of the bromodomain and extra-terminal domain (BET) family.<sup>1</sup> It is a ubiquitously expressed nuclear protein with roles in a variety of cellular processes, including regulation of gene transcription, cell cycle progression, and viral genome segregation.<sup>1,2</sup> BRD4 is comprised of two N-terminal bromodomains (BD1 and BD2) that can bind to acetylated lysine residues in histones, serving to couple histone acetylation marks to the transcriptional regulation of target promoters, an extra-terminal (ET) domain that facilitates protein-protein interactions, and a C-terminal motif, and can be expressed as either a long or short isoform generated *via* alternative splicing.<sup>1-3</sup> In addition to binding acetylated lysine residues on histones, BRD4 can bind to a variety of non-histone proteins and protein complexes, including positive transcription elongation factor b (P-TEFb) and 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>1,4-6</sup> In mice, knock-down of *Brd4* is embryonic lethal.<sup>2</sup> Chromosomal translocations leading to the fusion of *BRD4* with the nuclear protein in testis (*NUT*) gene and resulting in expression of BRD4-NUT fusion proteins are associated with NUT midline carcinoma (NMC).<sup>1,3</sup> Inhibition of BRD4-NUT binding to chromatin by the BET bromodomain inhibitor (+)-JQ1 (Item No. 11187) induces tumor regression and prolongs survival in NMC mouse xenograft models.<sup>7</sup> Cayman's BRD4 bromodomains 1 and 2 (human, recombinant; aa 2-477) 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. Wu, S.Y. and Chiang, C.M. The double bromodomain-containing chromatin adaptor Brd4 and transcriptional regulation. *J. Biol. Chem.* **282(18)**, 13141-13145 (2007).
3. Wang, C.Y. and Filippakopoulos, P. Beating the odds: BETs in disease. *Trends Biochem. Sci.* **40(8)**, 468-479 (2015).
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. Goldstein, D. P. and Kosasa, T. S. The subunit radioimmunoassay for human chorionic gonadotropin - clinical applications. *Progress in Gynecology* 145-184 (1975).

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