

PRODUCT INFORMATION



BRD3 bromodomains 1 and 2 (human, recombinant)

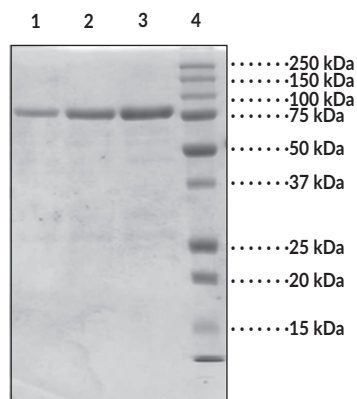
Item No. 14864

Overview and Properties

Synonyms:	Bromodomain containing protein 3, ORFX, RING3L, RING3-like protein
Source:	Recombinant N-terminal GST-tagged protein expressed in <i>E. coli</i>
Amino Acids:	2-434
Uniprot No.:	Q15059
Molecular Weight:	75.0 kDa
Storage:	-80°C (as supplied)
Stability:	≥2 years
Purity:	batch specific (≥85% estimated by SDS-PAGE)
Supplied in:	50 mM Tris, pH 8.0, with 150 mM sodium chloride and 20% glycerol
Protein Concentration:	batch specific mg/ml
Activity:	batch specific U/ml
Specific Activity:	batch specific U/mg

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

Image



Lane 1: BRD3 (4 µg)
Lane 2: BRD3 (6 µg)
Lane 3: BRD3 (8 µg)
Lane 4: MW Markers

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, 12/13/2021

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 protein 3 (BRD3) is a member of the bromodomain and extra-terminal domain (BET) family.¹ It 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. BRD3 is ubiquitously expressed and localizes to the nucleus where it preferentially binds acetylated histone H4 lysine 5 (H4K5), H4K12, and H4K14 at active transcription sites.^{2,3} Knockdown of *Brd3* enhances differentiation of C2C12 skeletal muscle myoblasts.⁴ Chromosomal translocations leading to the fusion of BRD3 with the nuclear protein in testis (NUT) gene and resulting in expression of BRD3-NUT fusion proteins are associated with NUT midline carcinoma (NMC).⁵ Cayman's BRD3 bromodomains 1 and 2 protein can be used for Western blot (WB).

References

1. Weidner-Glunde, M., Ottinger, M., and Schulz, T.F. WHAT do viruses BET on? *Front. Biosci.* **15**, 537-549 (2010).
2. Thorpe, K.L., Gorman, P., Thomas, C., et al. Chromosomal localization, gene structure and transcription pattern of the *ORFX* gene, a homologue of the MHC-linked *RING3* gene. *Gene* **200(1-2)**, 177-183 (1997).
3. LeRoy, G., Rickards, B., and Flint, S.J. The double bromodomain proteins Brd2 and Brd3 couple histone acetylation to transcription. *Mol. Cell* **30(1)**, 51-60 (2008).
4. Roberts, T.C., Etxaniz, U., Dall'Agnese, A., et al. BRD3 and BRD4 BET bromodomain proteins differentially regulate skeletal myogenesis. *Sci. Rep.* **7(1)**, 6153 (2017).
5. French, C.A., Ramirez, C.L., Kolmakova, J., et al. BRD-NUT oncoproteins: A family of closely related nuclear proteins that block epithelial differentiation and maintain the growth of carcinoma cells. *Oncogene* **327(15)**, 2237-2242 (2008).

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