

# PRODUCT INFORMATION



## CBX2 chromodomain (human recombinant)

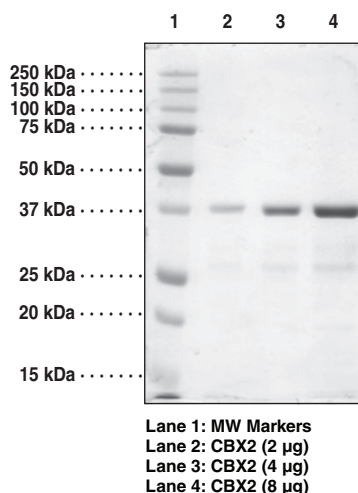
Item No. 14769

### Overview and Properties

**Synonyms:** CDCA6, Cell Division Cycle Associated 6, Chromobox Protein Homolog 2 (*Drosophila* Pc Class), M33, MGC10561, Modifier 3, SRXY5  
**Source:** Recombinant N-terminal GST-tagged protein expressed in *E. coli*  
**Amino Acids:** 2-90  
**Molecular Weight:** 38.2 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥6 months  
**Purity:** ≥90% estimated by SDS-PAGE  
**Supplied in:** 50 mM TRIS, pH 8.0, containing 150 mM sodium chloride and 20% glycerol  
**Protein Concentration:** *batch specific* mg/ml

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

### Image



**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, 06/21/2021

**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

---

CBX2 is a component of the polycomb repressive complex 1 and represses protein expression.<sup>1</sup> The chromodomain regions of CBX2 recognizes histone H3 trimethyl lysine 9 (H3K9me3) or trimethyl lysine 27 (H3K27me3).<sup>2-4</sup> CBX2 knock-out mice have impaired spleen formation, T cell expansion, and inactivation of the X chromosome, which can be lethal in some cases.<sup>5</sup> This protein product contains the chromodomain region of CBX2.

## References

---

1. Schoorlemmer, J., Marcos-Gutiérrez, C., Were, F., *et al.* Ring1A is a transcriptional repressor that interacts with the Polycomb-M33 protein and is expressed at rhombomere boundaries in the mouse hindbrain. *EMBO Journal* **16**(19), 5930-5942 (1997).
2. Nielsen, A.L., Oulad-Abdelghani, M., Ortiz, J.A., *et al.* Heterochromatin formation in mammalian cells: Interaction between histones and HP1 proteins. *Mol.Cell.* **7**(4), 729-739 (2001).
3. Bernstein, E., Duncan, E.M., Masui, O., *et al.* Mouse polycomb proteins bind differentially to methylated histone H3 and RNA and are enriched in facultative heterochromatin. *Mol. Cell Biol.* **26**(7), 2560-2569 (2006).
4. Kaustov, L., Ouyang, H., Amaya, M., *et al.* Recognition and specificity determinants of the human Cbx chromodomains. *J. Biol. Chem.* **286**(1), 521-529 (2011).
5. Katoh-Fukui, Y., Owaki, A., Toyama, Y., *et al.* Mouse polycomb M33 is required for splenic vascular and adrenal gland formation through regulating Ad4BP/SF1 expression. *Blood* **106**(5), 1612-1620 (2005).

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