

# PRODUCT INFORMATION



## JMJD2C tudor domain (human, recombinant)

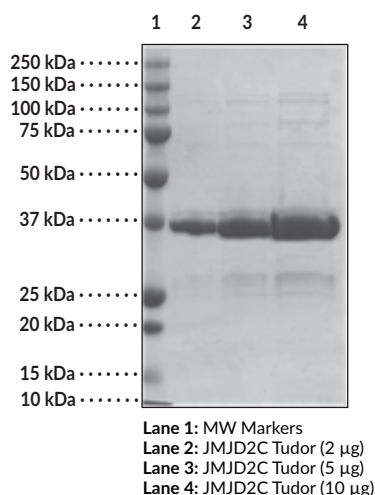
Item No. 9001953

### Overview and Properties

<b>Synonyms:</b>	GASC-1 Protein, KDM4C, Lysine-specific Demethylase 4C, JmjC Domain-containing Histone Demethylation Protein 3C, Jumonji Domain-containing Protein 2C
<b>Source:</b>	Recombinant N-terminal GST-tagged protein expressed in <i>E. coli</i>
<b>Amino Acids:</b>	869-1,003 (tudor domain)
<b>Uniprot No.:</b>	Q9H3R0
<b>M<sub>r</sub>:</b>	39.89 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥6 months
<b>Supplied in:</b>	50 mM Tris, pH 7.5, containing 150 mM sodium chloride and 20% 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



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, 06/18/2021

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

---

Tudor domains are small protein structural motifs of ~50 amino acids related to the “royal family” of methyl readers, which also includes chromo, MBT, PWWP, and Agenet-like domains.<sup>1,2</sup> Tudor domains occur either alone, in tandem, or with other domains and are found in many proteins that are involved in RNA metabolism, germ cell development, transposon silencing, DNA damage response, histone modification, and chromatin remodeling.<sup>3</sup> The tudor domains recognize symmetric methylated arginine or methylated lysine residues.<sup>4-7</sup> JMJD2C is an  $\alpha$ -ketoglutarate-dependent Fe (II) oxygenase that catalyzes the demethylation of trimethylated histone H3 at lysine residues 9 and 36 (H3K9me3 and H3K36me3).<sup>8</sup> This protein product contains the tandem tudor domains of JMJD2C.

## References

---

1. Maurer-Stroh, S., Dickens, N.J., Hughes-Davies, L., *et al.* The tudor domain ‘Royal Family’: Tudor, plant Agenet, Chromo, PWWP and MBT domains. *Trends Biochem. Sci.* **28(2)**, 69-74 (2003).
2. Lasko, P. Tudor domain. *Curr. Biol.* **20(16)**, R666-R667 (2010).
3. Chen, C., Nott, T.J., Jin, J., *et al.* Deciphering arginine methylation: Tudor tells the tale. *Nat. Rev. Mol. Cell Biol.* **12(10)**, 629-642 (2011).
4. Kim, J., Daniel, J., Espejo, A., *et al.* Tudor, MBT and chromo domains gauge the degree of lysine methylation. *EMBO Rep.* **7(4)**, 397-403 (2006).
5. Huang, Y., Fang, J., Bedford, M.T., *et al.* Recognition of histone H3 lysine-4 methylation by the double tudor domain of JMJD2A. *Science* **312(5774)**, 748-751 (2006).
6. Lee, J., Thompson, J.R., Botuyan, M.V., *et al.* Distinct binding modes specify the recognition of methylated histones H3K4 and H4K20 by JMJD2A-tudor. *Nat. Struct. Mol. Biol.* **15(1)**, 109-111 (2008).
7. Sprangers, R., Groves, M.R., Sinning, I., *et al.* High-resolution X-ray and NMR structures of the SMN tudor domain: Conformational variation in the binding site for symmetrically dimethylated arginine residues. *J. Mol. Biol.* **327(2)**, 507-520 (2003).
8. Pedersen, M.T., Agger, K., Laugesen, A., *et al.* The demethylase JMJD2C localizes to H3K4me3-positive transcription start sites and is dispensable for embryonic development. *Mol. Cell. Biol.* **34(6)**, 1031-1045 (2014).