

PRODUCT INFORMATION



TP53BP1 tudor-like region (human recombinant)

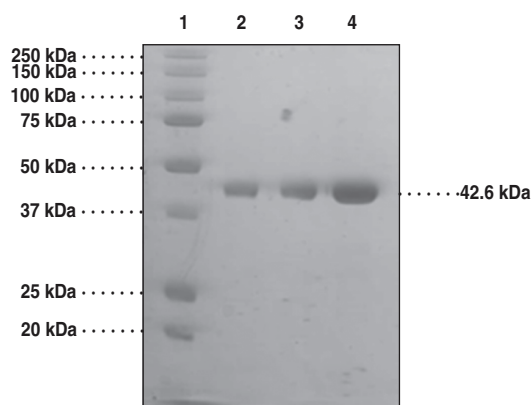
Item No. 14073

Overview and Properties

Synonyms: 53BP1, Tumor Suppressor p53-binding Protein 1
Source: Recombinant N-terminal GST-tagged protein expressed in *E. coli*
Amino Acids: 1717-1972 (N- and C-terminal truncations)
Uniprot No.: Q12888
Molecular Weight: 42.6 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 50 mM Tris-HCl, pH 8.0, containing 150 mM sodium chloride and 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: TP53BP1 (1 µg)
Lane 3: TP53BP1 (2 µg)
Lane 4: TP53BP1 (5 µg)

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.^{1,2} Tudor domains occur either alone, in tandem, or with other domains. They are found in many proteins that are involved in RNA metabolism, germ cell development, transposon silencing, DNA damage response, histone modification, and chromatin remodeling.³ Tudor domains recognize symmetric methylated arginine or methylated lysine residues.⁴⁻⁷ TP53BP1 binds both histone and non-histone substrates.^{4,8,9} Binding of TP53BP1 to dimethylated lysine 382 on p53 (p53 K382me2), as well as histone H4K20me2, through the tudor domain, has been shown to be important for TP53BP1 localization to DNA double strand breaks.^{8,10} This product contains the tudor domain of TP53BP1.

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**, 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. Kachirskaia, I., Shi, X., Yamaguchi, H., *et al.* Role for 53BP1 tudor domain recognition of p53 dimethylated at lysine 382 in DNA damage signaling. *J. Biol. Chem.* **283(50)**, 34660-34666 (2008).
9. Stewart, G.S. The demise of a TUDOR under stress opens a chromatin link to 53BP1. *EMBO J.* **31(8)**, 1847-1849 (2012).
10. Zgheib, O., Pataky, K., Brugger, J., *et al.* An oligomerized 53BP1 tudor domain suffices for recognition of DNA double-strand breaks. *Mol. Cell. Biol.* **29(4)**, 1050-1058 (2009).