

PRODUCT INFORMATION



WDR5 (human, recombinant)

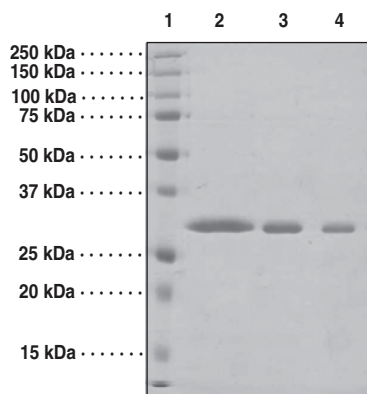
Item No. 10944

Overview and Properties

Synonyms:	BIG3, BMP2-induced 3-kb Gene Protein, SWD3, Set1c WD40 repeat protein, homolog, WD-Repeat Protein 5
Source:	Recombinant protein expressed in <i>E. coli</i> . An N-terminal hexahistidine tag and SUMOpro tag were removed by cleavage with SUMO Protease 1 (Ulp1). <i>SUMOpro and SUMO Protease 1 were used under non-exclusive license from LifeSensors, Inc. www.lifesensors.com.</i>
Amino Acids:	2-334 (full length)
Molecular Weight:	34.4 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability:	≥1 year
Purity:	<i>batch specific</i> (≥95% estimated by SDS-PAGE)
Supplied in:	50 mM Tris, pH 8.0, with 150 mM sodium chloride and 20% glycerol
Protein	
Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg

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

Image



Lane 1: MW Markers
Lane 2: WDR5 (4 µg)
Lane 3: WDR5 (2 µg)
Lane 4: WDR5 (1 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, 02/28/2020

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

WD repeats are motifs of approximately 40 amino acids typically terminating in conserved tryptophan and aspartate residues. Several WD40 repeats combine to form the structural WD domain. Such an arrangement facilitates protein-protein interactions allowing the formation of multiprotein complexes.¹ Human WDR5 contains seven WD 40 repeats and was originally identified to accelerate osteoblast differentiation.² WDR5 has been demonstrated to bind histone H3 by recognizing the first three amino acids of the N-terminal tail.³ Binding of WDR5 to a conserved arginine-containing motif in MLL-1, the so-called WDR5 interaction ("Win") motif, promotes the assembly and activity of the MLL core complex.⁴ Additional interactions have been demonstrated with other components of the human MLL core protein complex, which includes ASH2L and RbBP5.^{5,6} MLL1-5 protein complexes catalyze the di- and trimethylation of histone H3 at lysine 4 (H3K4me2/me3), leading to the maintenance of global H3K4 trimethylation.⁷

References

1. Neer, E.J., Schmidt, C.J., Nambudripad, R., *et al.* The ancient regulatory-protein family of WD-repeat proteins. *Nature* **371(6495)**, 297-300 (1994).
2. Gori, G., Divieti, P., and Demay, M.B. Cloning and characterization of a novel WD-40 repeat protein that dramatically accelerates osteoblastic differentiation. *J. Biol. Chem.* **276(49)**, 46515-22 (2001).
3. Couture, J.-F., Collazo, E., and Trievel, R.C. Molecular recognition of histone H3 by the WD40 protein WDR5. *Nat. Struct. Mol. Biol.* **13(8)**, 698-703 (2006).
4. Patel, A., Vought, V.E., Dharmarajan, V., *et al.* A conserved arginine-containing motif crucial for the assembly and enzymatic activity of the mixed lineage leukemia protein-1 core complex. *J. Biol. Chem.* **283(47)**, 32162-75 (2008).
5. Odho, Z., Southall, S.M., and Wilson, J.R. Characterization of a novel WDR5-binding site that recruits RbBP5 through a conserved motif to enhance methylation of histone H3 lysine 4 by mixed lineage leukemia protein-1. *J. Biol. Chem.* **285(43)**, 32967-76 (2010).
6. Dou, Y., Milne, T.A., Ruthenburg, A.J., *et al.* Regulation of MLL1 H3K4 methyltransferase activity by its core components. *Nat. Struct. Mol. Biol.* **13(8)**, 713-9 (2006).
7. Wang, P., Lin, C., Smith, E.R., *et al.* Global analysis of H3K4 methylation defines MLL family member targets and points to a role for MLL1-mediated H3K4 methylation in the regulation of transcriptional initiation by RNA polymerase II. *Mol. Cell. Biol.* **29(22)**, 6074-85 (2009).

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