

Data Sheet

MLL1/WDR5/Ash2L/ RbBP5 Complex

Human, recombinant

Catalog #: 51022

Lot #: 120906 **Conc.:** 0.24 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.4, 300 mM NaCl, 250 mM imidazole, 1 mM TCEP, 1 μ M ZnCl₂, and 10% glycerol.

Stability: >6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

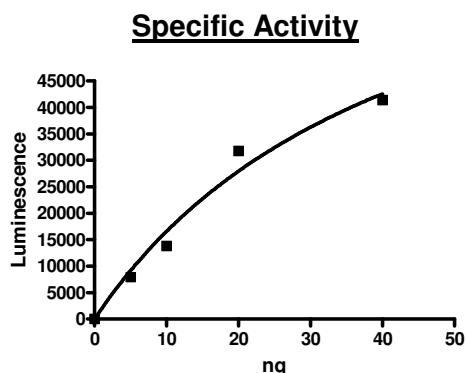
1. Patel, A., *et al.*, *J. Biol. Chem.* 2009; **284(36)**: 24242-24256.
2. Mohan, M., *et al.*, *Nat. Rev. Cancer* 2010; **10(10)**: 721-728.

Description: MLL1/WAR complex of human **MLL1** (GenBank Accession No. NM_005933) a.a. 3745-end with N-terminal His-tag and MW= 27 kDa, human **Ash2L** (GenBank Accession No. NM_001105214), a.a. 2-end with N-terminal His-tag and MW= 61 kDa, and human **RbBP5** (GenBank Accession No. NM_005057), a.a. 2-end with N-terminal His-tag and MW= 60 kDa individually expressed in an *E. coli* expression system, and human **WDR5** (GenBank Accession No. NM_017588), a.a. 2-end with N-terminal His-tag and MW= 37.4 kDa expressed in a baculovirus Sf9 expression system.

Assay Conditions: 50 μ l reaction mix (50 mM Tris, pH 8.8, 5 mM MgCl₂, 4 M DTT, 20 μ M S-adenosylmethionine, and 5-50 ng MLL complex) add to the wells coated with the substrate. Incubate at room temperature for 1 hr. Add antibody against methylated K4 residue of histone H3, incubate 1 hr. Then, incubate 30 minutes with the secondary HRP-labeled antibody. Finally, add HRP chemiluminescent substrate and read luminescence.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**4-20% SDS-PAGE
Coomassie Staining**

Lane 1:

MLL1/WAR Complex

Lane 2:

Protein Marker

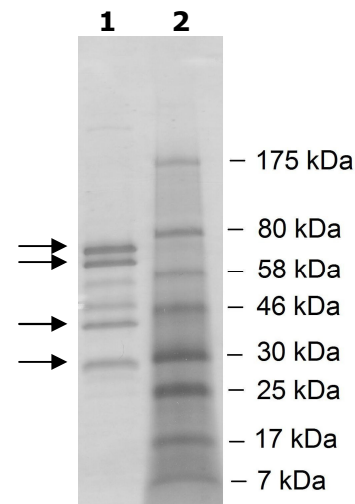
Purity: \geq 82%

MW: RbBP5: 60 kDa

Ash2L: 61 kDa

WDR5: 37 kDa

MLL1: 27 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Sequences:

MLL1

MHHHHHHFRFHKPEEANEPLNPHGSARAEVHLRKSADFDMFNFLASKHRQPPEYNP
NDEEEEEEVQLKSARRATSMPLPMPFRFRHLKKTKEAVGVYRSPHGRGLFCKRND
AGEMVIEYAGNVIRSIQTDKREKYYDSKGIGCYMFRIDDSEVVDATMHGNAARFINHS
CEPNCYSRVINIDGQKHIVIFAMRKIYRGEELTYDYKFPIEDASNKLPCNCGAKKCRKF
LN

WDR5

MHHHHHHAATEKKPETEAARAQTPSSSATQSKPTPVKPNYALKFTLAGHTKAVSSV
KFSNGEWLASSSADKLIKWGAYDGKFEKTISGHKLGISDVAWSSDSNLLVSASDDK
TLKIWDVSSGKCLKTLKGHSNYVFCCNFNPQSNLIVSGSFDESVRWDVKTGKCLKTL
PAHSDPVS AVHFNRDGLIVSSSYDGLCRIWDTASGQCLKTLIDDDNPPVSFVKFSPN
GKYILAAATLDNTLKLWDYSKGKCLKTYTGHNKCYCIFANFSVTGGKWIVSGSEDNLV
YIWNLQTKIVQKLQGHDTVVISTACHPTENIIASAALENDKTIKLWKSDC

Ash2L

MHHHHHHDTQAGSVDEENGRQLGEVELQCGICTKWFTADTFGIDTSSCLPFMTNYSF
HCNVCHHSGNTYFLRKQANLKEMCLSALANLTWQSRTQDEHPKTMFSKDKDIIPFID
KYWECMTTRQRP GKMTWPNNIVKTMSKERDVFLVKEHPDPGSKDPEEDYPKFGLLD
QDLSNIGPAYDNQKQSSAVSTSGNLNGGIAAGSSGKGRGAKRKQQDGGTTGTTKKA
RSDPLFSAQRLPPHGYPLEHPFNKDG YRYLAEPDPHAPDPEKLELDCWAGKPIPGD
LYRACL YERVLLALHDRAPQLKISDDRLTVVGEKGYSMVRASHGVRKGAWYFEITVD
EMPPDTAARLGWSQPLGNLQAPLG YDKFSYSWRSKKGTKFHQSIGKHYSSGYGGG
DVLGFYINLPEDTETAKSLPDTYKDKALIKFKSYLYFEEKDFVDKAEKSLKQTPHSEIIF
YKNGVNQGVAYKDIFEGVYFPAISLYKSC TVSINFGPCFKYPPKDLTYRPMSDMGWG
AVVEHTLADVLYHVETEVDGRRSPPEP

RbBP5

MHHHHHHNLELLESFGQNYPEEADGTLDCISMAL TCTFN RWG TLLAVGCNDGRIVW
DFLTRGIAKII SAHIHPVCSLCWSRDGHKLV SASTDNIVSQWDVLSGDCDQRFRFPSP
LKVQYHPRDQNKVLVCPMKSAPVMLT LSDSKHV VLPVDDSDLN VVASFDRRGEYIY
TGNAGKILVLKTDSDQLVASFRVTTGTSNTTAIKSIEFARKGSCFLINTADRIIRVYDG
REILTCGRDGEPEPMQKLQDLVNRTPWKKCCFSGDGEYIVAGSARQHALYIWEKSIG
NLVKILHGTRGELLLDVAWHPVRPIIASISSGVVSIWAQNQVENWSAFAPDFKELDEN
VEYEERESEFDIEDEDKSEPEQTGADAAEDEVDVTSVDPIAAFCSSEDELEDKALL
YLP I APEVEDPEENPYGPPDAVQTS LMDEGASSEK KRQSSADGSQPPKKPKTTNI
ELQGV PNDEVHPLLGVKGDGKSKKKQAGR PKGSKGKEKDS PFKPKLYKGDRGLPL
EGSAKGVQAELSQPLTAGGAISELL

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com