

## Data Sheet

### RbBP5 (2-end)

Human, recombinant, N-terminal His tag

**Catalog #:** 71202

**Lot #:** 120206-G      **Conc.:** 0.22 mg/ml

**Formulated in:** 20 mM Tris-HCl, pH 7.4,  
300 mM NaCl, 10% Glycerol, 1 mM TCEP,  
and 1  $\mu$ M ZnCl<sub>2</sub>.

**Stability:** >6 months at -80°C

**References:**

1. Cao, F., *et al.*, *PLoS One*. 2012 Nov 23; **5(11)**: e14102.
2. Odho, Z., *et al.*, *J Biol Chem*. 2010 Oct 22; **285(43)**: 32967-76.

**Description:**

Human retinoblastoma binding protein 5 (RbBP5), GenBank Accession No. NM\_005057, a.a. 2 – 538 (end) with N-terminal His-tag, MW= 60 kDa, expressed in an *E. coli* expression system.

**Application:**

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

## Quality Assurance

**10% SDS-PAGE  
Coomassie  
staining**

**Lane 1:**

4  $\mu$ g RbBP5

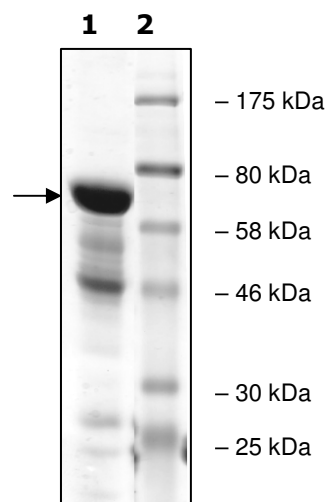
**Lane 2:**

Protein Marker

BioLabs (#P7708L)

**MW:** 60 kDa

**Purity:**  $\geq$ 65%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

**Protein Sequence:**

**MHHHHH**NLELLESFGQNYPEEADGTLDCISMALCTFNRWGTL LAVGCNDGRIVW  
DFLTRGIAKIISAHHPVCSLCWSRDGHKLV SASTDNIVSQWDVLSGDCDQRF RFPSP  
LKVQYHPRDQNKVLVCPMKSAPVMLT LSDKHVLPVDDSDLN VVASFDRRGEYIY  
TGNAGKILVLKTDSQDLVASFRVTTGTSNTT AIKSIEFARKGSCFLINTADRIIRVYDG  
REILTCGRDGEPEPMQKLQDLVNRTPWK KCCFSGDGEYIVAGSARQHALYIWEKSIG  
NLVKILHGTRGELL LDVAWHPVRPIIASISSGVVSIWAQNQVENWSAFAPDFKELDEN  
VEYEERESEFDIEDEKSEPEQTGADAAE DEEVDVTSVDPIAAFCSSEDELEDKALL  
YLP IAPEVEDPEENPYGPPDAVQTSLMDE GASSEKQRQSSADGSQPPKKPKTTNI  
ELQGVPNDEVHPLLGVKGDGKSKKKQAGR PKGSKGKEKDSPFKPKLYKGDRGLPL  
EGSAKGKVQAELSQPLTAGGAISELL

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)