

## Data Sheet

### PHIP (1146-1287)

Human, recombinant, N-terminal His-tag

**Catalog #:** 31140

**Lot #:** 130822      **Conc.:** 0.65 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Stability:** >6 months at  $-70^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

1. De Semir, D., *et al.*, *Proc Natl Acad Sci USA*. 2012 May 1; **109(18)**:7067-72.
2. Farhang-Fallah, J., *et al.*, *J Biol Chem*. 2000 Dec 22; **275(51)**:40492-7.

**Description:**

Human bromodomain pleckstrin homology domain interacting protein (PHIP), also known as IRS-1 PH Domain-Binding Protein, WDR11, and BRWD2, GenBank Accession No. NM\_017934, a.a. 1146 - 1287, with N-terminal His-tag, MW = 17 kDa, expressed in an *E. coli* expression system.

**Application:**

Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

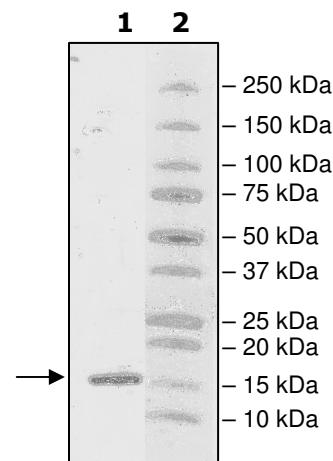
## Quality Assurance

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

**Lane 1:**  
2.3  $\mu\text{g}$  PHIP

**Lane 2:**  
Protein Marker

**MW:** 17 kDa  
**Purity:**  $\geq 90\%$



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 West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

MHHHHHSLIYKPLDGEWGTNPRDEECERIVAGINQLMTLDIASAFVAPVDLQAYPMY  
CTVVAYPTDLSTIKQRLNRFYRRVSSLMWEVRYIEHNTRTFNEPGSPIVKSASFVTDL  
LLHFIKDQTCYNIIPLYNSMKKKVLSDSEDEE

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)