

Data Sheet

PHIP (1146-1287)

Human, recombinant, N-terminal GST-tag

Catalog #: 31142

Lot #: 130724-G1 Conc.: 0.10 mg/ml

Formulated in: 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol.

Stability: >6 months at -70°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 GST-tag, MW = 43 kDa, expressed in an *E. coli* expression system.

Application:

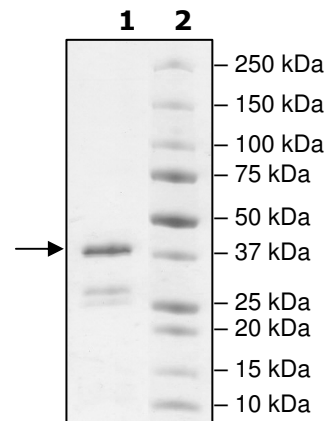
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg PHIP
Lane 2:
Protein Marker

MW: 43 kDa
Purity: $\geq 71\%$



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 West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPQIDKYLKSSKYIAWPLQGQATFGGGDHPPKSD**LEVLFGGPLGSSLIYK**
PLDGEWGTNPRDEECERIVAGINQLMTLDIASAFVAPVDLQAYPMYCTVVAYPTDLST
IKQRLNRFYRRVSSLMWEVRYIEHNTRTFNEPGSPIVKS AKFVTDLLLHFIKDQTCYNI
IPLYNSMKKKVLSDEDEE

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