

Data Sheet

WDR9 (1308 – 1436)

Human, recombinant, N-terminal GST-tag

Catalog #: 31135

Lot #: 130719-D Conc.: 3.75 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, and 20% glycerol.

Stability: >6 months at –80°C. **Avoid
multiple freeze/thaw cycles.**

References:

1. Huang, H., *et al.*, *Dev Dyn.* 2003 Aug;
227(4): 608-14.
2. Ramos, VC., *et al.*, *Biochim Biophys
Acta.* 2002 Sep 27;**1577(3):** 377-83.

Description:

Human bromodomain and WD repeat domain containing 1 (BRWD1) or WDR9, amino-acids 1308 - 1436, (GenBank Accession No. NM_018963) with N-terminal GST-tag, MW = 42 kDa, expressed in an *E. coli* expression system.

Application:

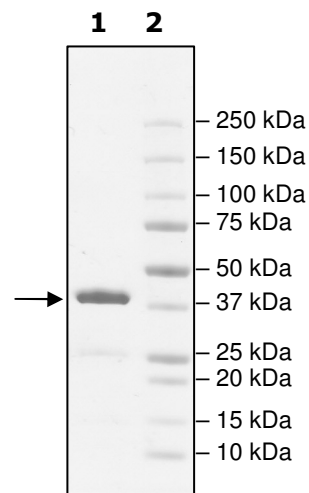
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3 µg WDR9
Lane 2:
Protein Marker

MW: 42 kDa
Purity: ≥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
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:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPQIDKYLKSSKYIAWPLQGQATFGGGDHPPKSDLEVLFGGPLGSIRATN
YVESNWKKQCKELVNLIFQCEDESEPFRRQPVDLVEYPDYRDIIDTPMDFGTVRETL DAG
NYDSPLEFCKDIRLIFSNAKAYTPNKRSKIYSMTLRLSALFEEKMKKISSDFKIGQKFNE
KLRRSQ

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