

## Data Sheet

### WDR5

Human, recombinant, N-terminal His-tag

**Catalog #:** 71200

**Lot#:** 120912      **Conc.:** 0.50 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 200 mM  
Imidazole, and 20% glycerol.

**Stability:** >6 months at -80°C

#### References:

1. Trievel, R.C., Shilatifard, A. (2009). *Nature Struct. Mol. Biol.* **16 (7)**: 678-80.
2. Wang, Y.Y., et al. (2010). *Proc. Natl. Acad. Sci. USA* **107(2)**:815-20.
3. Odho, Z., et al. (2010). *J. Biol. Chem.* **285(43)**:32967-76.

#### Description:

Human WDR5 (WD Repeat Domain 5), also known as SWD3 or BIG-3, GenBank Accession No. NM\_017588, a.a. 2-334(end), with N-terminal His-tag, MW=37.5 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

#### Application:

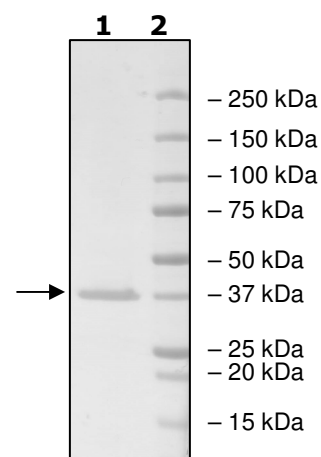
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

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

**Lane 1:**  
2 µg WDR5  
**Lane 2:**  
Protein Marker

**MW:** 38 kDa  
**Purity:** ≥95%



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)