

Data Sheet

WDR5, His-Tag, Biotin-Labeled

Human, Recombinant, N-terminal Avi-His-tag

Catalog #: 80016

Lot #: 160902

Conc.: 0.16 mg/ml

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

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended. If necessary, use carrier protein (BSA 0.1-0.5%).*

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 Avi-His-tag, MW=39 kDa, expressed in Sf9 insect cells via a baculovirus expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$.

Application:

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

Quality Assurance

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

Lane 1:

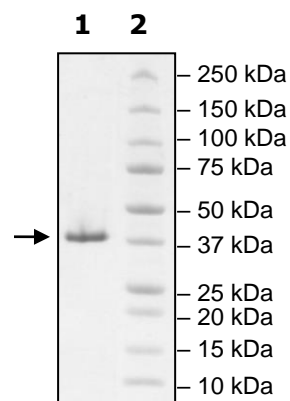
WDR5

Lane 2:

Protein Marker

MW: 39 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

Please visit our website at: www.bpsbioscience.com