

Data Sheet

DDX58, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100070

Lot #: 180723

Conc.: 0.69 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3mM DTT, 100 µg/ml flag peptide

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. Jang, M., *et al.*, *Cell*. 2015; **96(2)**: 266-274.
2. Suarez-Calvet, X., *et al.*, *Journal of Pathology*. 2014; **233(3)**: 258-268.

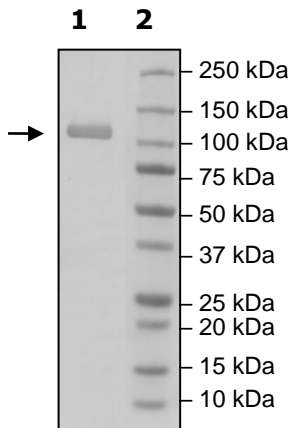
Description: Human DDX58 protein, also known as DEAD box protein 58, retinoic acid inducible gene 1, and RIG-1, GenBank Accession No. NM_014314, a.a. 2-925 (end), with N-terminal FLAG-tag, MW=108 kDa, expressed in a HEK293 expression system.

Application: Useful as a positive control for Western blot studies.

Quality Assurance

**Western Blot
w/ Flag-antibody**

Lane 1:
DDX58
Lane 2:
Protein Marker

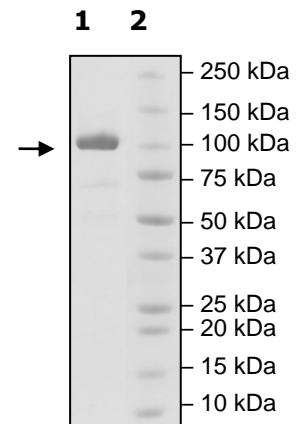


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
DDX58
Lane 2:
Protein Marker

Purity: ≥95%

MW: 108 kDa



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