

## Product Information

|                           |  |
|---------------------------|--|
| <b>Construct:</b>         | DDX58 (FLAG-2-925(end))  |
| <b>Concentration:</b>     | 0.73 mg/ml   |
| <b>Species:</b>           | Human  |
| <b>Formulated In:</b>     | 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 µg/ml FLAG peptide, 20% glycerol, 3 mM DTT |
| <b>Expression System:</b> | HEK293   |
| <b>Format:</b>            | Aqueous buffer solution  |
| <b>Stability:</b>         | At least 6 months at -80°C. Avoid freeze/thaw cycles.  |
| <b>Storage:</b>           | -80°C  |
| <b>Genbank Accession:</b> | NM_014314  |
| <b>MW:</b>                | 108 kDa  |
| <b>Purity:</b>            | ≥90%   |

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining

