

Product Information

| | |
|---------------------------|---|
| Construct: | HSD17B13 (1-300(end)-FLAG) |
| Concentration: | 0.13 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 100 µg/ml FLAG peptide |
| Expression System: | HEK293 |
| Format: | Aqueous buffer solution |
| Stability: | At least 6 months at -80°C. Avoid freeze/thaw cycles. |
| Storage: | -80°C |
| Genbank Accession: | NM_178135 |
| MW: | 37 kDa |
| Purity: | 86% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

