

Product Information

| | |
|---------------------------|--|
| Construct: | NLRP3 (FLAG-Pre-134-1036(end)) / NEK7 (His-Pre-2-302(end)) |
| Concentration: | 0.33 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 µg/ml FLAG peptide, 3 mM DTT, and 20% glycerol |
| Expression System: | HEK293 |
| Format: | Aqueous buffer solution |
| Stability: | At least 6 months at -80°C. Avoid freeze/thaw cycles. |
| Storage: | -80°C |
| Genbank Accession: | NLRP3: NM_004895; NEK7: NM_133494 |
| MW: | NLRP3: 105 kDa; NEK7: 37 kDa |
| Purity: | ≥90% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

