

Product Information

Construct:	NLRP3 (delPYD) (FLAG-134-1036(end))
Mutation:	delPYD
Concentration:	0.17 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 500 mM KCl, 0.5 mM TCEP, and 10% 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:	NM_004895
MW:	105 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

