

NLRP3 (NALP3), His-FLAG-Tags (Sf9-derived) Recombinant

Catalog: 80117
Lot: 230323-A

Product Information

Description:	Human NLR Family Pyrin Domain Containing 3 (NLRP3, also known as NALP3), full length encompassing amino acids 2-1036(end). This construct contains an N-terminal His-tag (6xHis) followed by a-FLAG-tag. The recombinant protein was affinity purified.
Species	Human
Construct:	NLRP3 (His-FLAG-2-1036(end))
Concentration:	1.24 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 500 mM KCl, 0.5 mM TCEP, 20% glycerol, and 100 µg/ml FLAG peptide.
MW:	120 kDa
Genbank Accession:	NM_004895
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

