

# NLRP3 (K232A), His-FLAG-Tag (Sf9-derived) Recombinant

Catalog: 40741  
Lot: 221027

## Product Information

<b>Construct:</b>	NALP3 (K232A) (His-FLAG-2-1036(end))
<b>Mutation:</b>	K232A
<b>Concentration:</b>	0.23 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 µg/ml FLAG peptide, and 20% glycerol.
<b>Expression System:</b>	Sf9
<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_004895
<b>MW:</b>	120 kDa
<b>Purity:</b>	58%

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

