

## Data Sheet

### NLRP3 (NALP3), His-FLAG-Tags

Human, Recombinant, N-terminal  
His-FLAG-tags

**Catalog #:** 100189

**Lot #:** 190426-2

**Conc.:** 0.14 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 0.8 M Arginine, 100 µg/ml FLAG peptide, 3 mM DTT and 20% glycerol.

**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles.

**Description:** Human NLR Family Pyrin Domain Containing 3 (NLRP3), also known as NALP3, Cold Autoinflammatory Syndrome 1 Protein (CIAS1), cryopyrin and PYPAF1, Genbank Accession No. NM\_004895, a.a. 2-1036(end) with N-terminal His-FLAG-tag, expressed in a HEK293 expression system. MW = 120 kDa.

#### References:

1. Kummer, J.A., *J. Histochem. Cytochem.*, 2007 May; **55(5)**:443-52.
2. Manji, G.A., *J. Biol. Chem.* 2002 Mar 29; **277(13)**:11570-5.

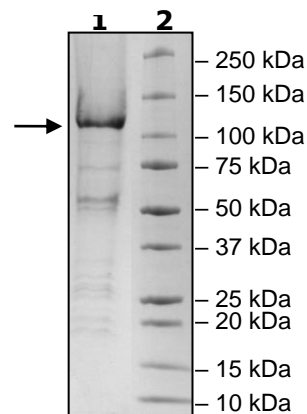
## Quality Assurance

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

**Lane 1:**  
2µg NLRP3  
**Lane 2:**  
Protein Marker

**MW:** 120 kDa

**Purity:** ≥75%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)