

Data Sheet

NLRP3 (NALP3) (GKT(231-233)AAA), His-FLAG-Tag

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

Catalog #: 100190

Lot #: 190423

Conc.: 0.28 mg/ml

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, 3mM DTT and 20% glycerol.

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

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.

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 and with GKT→AAA mutations at 231-233, expressed in a HEK293 expression system. MW = 120 kDa.

Quality Assurance

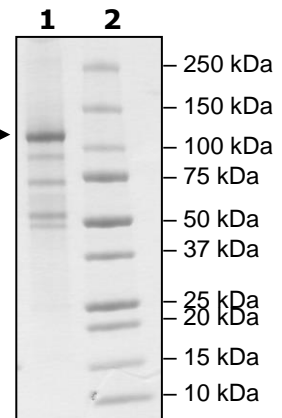
4-20% SDS-PAGE Coomassie staining

Lane 1:
2µg NLRP3

Lane 2:
Protein Marker

MW: 120 kDa

Purity: ≥50%



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

Please visit our website at: www.bpsbioscience.com