Catalog # **Aliquot Size**

20 µg

100 µg

C19NP-60DH-20 C19NP-60DH-100

Anti 2019-nCoV N-Protein hlgG Antibody, HRP conjugated

Recombinant Human Monoclonal IgG Antibody

Cataloa # C19NP-60DH

Lot # X3397-14

Cited Applications

ELISA, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the 2019-nCoV N-Protein

Cross Reactivity

SARS-CoV N-Protein

Host/Isotype/Clone

Human laG

Formulation

1X PBS + 50% alycerol + HRP stabilizer

Alternative Name(s)

Anti-SARS-CoV 2 N-Protein antibody

Stability

1yr at -20°C from date of shipment

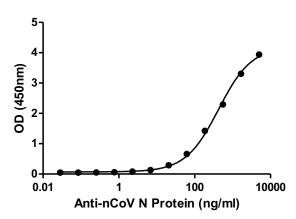
Scientific Background

nCoV N-Protein / Nucleocapsid protein is an RNA-binding protein primarily involved in binding to the viral RNA genome during packaging of viral RNA into a helical ribonucleoprotein or nucleocapsid structure (1). In addition, N-protein regulates viral RNA synthesis during replication/transcription and modulates host-pathogen interactions, such as actin reorganization, host cell cycle progression, and apoptosis (2). N-Protein is abundantly expressed during infection and antibodies against N-Protein have been detected in confirmed cases of Covid-19. Therefore, N-Protein is a potent antigen that provides basis for future vaccine and diagnostic kit development (3).

References

- McBride R, et al: The coronavirus nucleocapsid is a multifunctional protein. Viruses. 2014, 6:2991-3018.
- Surjit M, et al: The nucleocapsid protein of severe acute respiratory syndrome-coronavirus inhibits the activity of cyclin-cyclin-dependent kinase complex and blocks S phase progression in mammalian cells. J Biol Chem. 2006, 281:10669-10681.
- Zeng W, et al: Biochemical characterization of SARS-CoV-2 nucleocapsid protein. Biochem Biophys Res Commun. 2020, 527(3):618-623.

Assay Data



Binding of Anti 2019-nCoV N Protein hlgG Antibody, HRP conjugated (C19NP-60DH) to immobilized 2019-nCoV RBD-N Protein Chimera (C19NSD-G241H) determined by ELISA.

Anti 2019-nCoV N-Protein hlgG Antibody, **HRP** conjugated

Recombinant Human Monoclonal IgG Antibody

Cataloa# C19NP-60DH Lot # X3397-14 Concentration 0.2 μg/μl

Stability

1yr at -20°C from date of shipment Store product at -20°C. For optimal Storage & Shipping

storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com or IVD@signalchem.com - www.signalchem.com

Revised date: 2022-09-07 Page 1 of 3

SAFETY DATA SHEET

Article 1 - Product Identification

Product Name: Anti 2019-nCoV N-Protein hlgG Antibody, HRP conjugated Catalog # C19NP-60DH

This product is sold only for research use by qualified laboratory personnel, and is not to be used as a drug, medical device, food additive, cosmetic, nor household chemical. It is not to be used in diagnostic, therapeutic, consumer, agricultural, nor pesticidal applications.

Manufacturer's Name: SignalChem Biotech Inc. Street Address: 110-13120 Vanier Place City, Prov. Postal Code: Richmond, BC, V6V 2J2

Fax: 604-232-4601 EMERGENCY PHONE: 604-232-4600

Article 2 - Hazard Identification

WHMIS/GHS classification: This mixture is self-classified as non-hazardous. It does not contain hazardous material at a
concentration that would be considered hazardous.

Signal words: None.
 Hazard statements: None.
 Precautionary statements:

P260: Do not breath dust or mist

P264: Wash hands thoroughly after handling.

P280: Wear protective gloves

P301+P330+P331: IF SWALLOWED: rinse mouth. Do Not induce vomiting.

P303+P361+P353: IF ON SKIN: Take off contaminated clothing. Rinse skin with water/shower.

P363: Wash contaminated clothing before reuse.

P304+P340: IF INHALED: Remove victim to fresh air and keep at rest in a position comfortable for breathing. P305+P351+P338: IF IN EYES: rinse cautiously with water for several minutes. Remove contact lenses if present

and safe to do so. Seek medical attention.

Article 3 - Composition/Information on Ingredients

Chemical Characterization: Mixtures.

Description: This product consists of the substances listed below.

Common name	Chemical name	CAS-No.	Concentration
Sodium Chloride	Sodium Chloride	7647-14-5	0.72%
Sodium Phosphate, Dibasic	Sodium Phosphate, Dibasic	7782-85-6	0.248%
2-Chloroacetamide	2-Chloroacetamide	79-07-2	<0.1%
Potassium Phosphate, Monobasic	Potassium Phosphate, Monobasic	7778-77-0	0.024%
Potassium Chloride	Potassium Chloride	7447-40-7	0.02%
Protein	-	-	0.02%

Article 4 - First-aid Measures

- General information: Consult a physician by providing the SDS.
- After inhalation: Move to fresh air. If cannot breathe, give artificial respiration and consult a physician.
- After skin contact: Remove contaminated clothing. Immediately wash with soap and plenty of water and rinse thoroughly. Wash contaminated clothing before re-use.
- After eye contact: Rinse opened eyes with plenty of water for at least 15 minutes. Remove contact lenses, if present and easy to do so. Consult a physician.
- After swallowing: If the patient is conscious, rinse the mouth with plenty of water and consult a physician. Do not induce vomiting. Never give anything by mouth to an unconscious person.

Revised date: 2022-09-07 Page 2 of 3

SAFETY DATA SHEET

Article 5 - Fire-fighting Measures

- Suitable extinguishing media: Use water spray, extinguishing powder, carbon dioxide, or other appropriate measure that is suitable to the environment.
- Specific hazards arising from the substance or mixture: None known.
- Special protective equipment and precautions for fire-fighters: Self-contained breathing apparatus if necessary.

Article 6 - Accidental Release Measures

- Personal precautions, protective equipment and emergency procedures: Apply standard laboratory practices and personal protective equipment. Avoid breathing vapors, mist, or gas. Ensure adequate ventilation.
- Environmental precautions: Do not allow to enter drains.
- Methods and materials for containment and cleaning up: Absorb on sand or vermiculite and place in closed containers for disposal.

Article 7 - Handling and Storage

- Precautions for sate handling: Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact
 with eyes, skin or clothing.
- Conditions for safe storage: Store in a dry and well-ventilated place in -20 °C. Keep container upright and tightly closed.

Article 8 - Exposure Controls/Personal Protection

- Components with limit monitoring values at workplace: N/A
- Appropriate engineering controls:
 - Apply adequate ventilation including mechanical exhaust or laboratory fume hood. Follow standard laboratory practices.
- Individual protection measures:
- Respiratory protection:
 - Use appropriate respirator if there is inadequate ventilation by following the government standards.
- Hand protection:
 - Wear gloves and use proper glove removal technique to avoid skin contact. Discard gloves after use by following the applicable laboratory regulations. Wash and dry hands.
- Eye/face protection:
 - Safety goggles with side-shields approved under appropriate government standards.
- Skin/body protection:
 - Use appropriate clothing, footwear and any additional protection measures to protect from splashing or contamination.

Article 9 - Physical and Chemical Properties

Appearance: Colorless fluid.	Danger of explosion: Product does not present an explosion hazard.	
Odour/Odour Threshold: Not determined.	Explosion limits: Not available.	
pH: Not available.	Decomposition temperature: Not available.	
Melting point/freezing point: Not determined.	Vapor pressure at 20 °C: Not determined.	
Boiling point/Boiling range: Not determined.	Density: Not determined.	
Flash point: Not determined.	Relative density: Not determined.	
Flammability (solid, gaseous): Not determined.	Vapor density: Not determined.	
Ignition temperature: Not determined	Evaporation rate: Not determined.	
Auto-igniting: Product is not self-igniting.	Solubility in / Miscibility with Water: Fully miscible.	

Article 10 - Stability and Reactivity

- Reactivity: Stable under recommended transport and storage conditions.
- Chemical stability: Stable under recommended transport and storage conditions.
- Possible hazardous reactions: No dangerous reactions known.
- Conditions to avoid: None determined.
- Incompatible materials: None known.
- Hazardous decomposition products: No hazardous decomposition products or polymerization under normal conditions.

Revised date: 2022-09-07 Page 3 of 3

SAFETY DATA SHEET

Article 11 - Toxicological Information

- Acute toxicity: Not available.
- LD/LC50: Sodium azide: None expected.
- Skin corrosion/irritation: None expected.
- Serious eye damage/eye irritation: None expected.
- Respiratory or skin sensitization: Not available.
- Germ cell mutagenicity: Not available.
- Carcinogenicity: No components are listed in IARC, or NTP, or OSHA, or ACGIH.
- Reproductive toxicity: Not available.
- Teratogenicity: Not available.
- Specific target organ toxicity single exposure/ repeated exposure (GHS): Not available.
- Aspiration hazard: Not available.
- Potential health effects:
 Inhalation: None expected.
 Ingestion: None expected.
 Skin: None expected.
 - Eyes: None expected.
- Signs and Symptoms of Exposure: Not available.
- Synergistic effects: Not available.

Article 12 - Ecological Information

- Eco-toxicity: Not applicable.
- Biodegradability: Not applicable.
- Bio-accumulative potential: Not applicable.
- Mobility in soil: Not applicable.
- PBT and vPvB assessment: Not applicable.
- Other adverse effects: Not applicable.

Article 13 - Disposal Considerations

- Disposal methods: In accordance to applicable national, regional, or local laws and regulations. For additional handling information and protection of employees please refer to Article 7 and 8.
- Contaminated packaging: Disposal should be made in accordance to official regulations. Use water or cleansing agents to clean the area.

Article 14 - Transport Information

- **DOT**: Not dangerous goods.
- IMDG: Not dangerous goods.
- IATA: Not dangerous goods.

Article 15 - Regulatory Information

- WHMIS Classification: Non-hazardous.
- GHS label elements: Not applicable.
- Signal word: Not applicable.
- Hazard statements: Not applicable.

Article 16 - Other Information

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. SignalChem shall not be held liable for any damage resulting from handling or from contact with the above product. See the Technical Specification, Packing Slip, Invoice, and Product Catalog for additional terms and conditions of sale.