

# ACE2 (19-740) Protein

Recombinant human protein expressed in CHO cells

#### Catalog # A51C2-G341F

Lot # O3814-4

# **Product Description**

Recombinant human angiotensin-converting enzyme 2 (ACE2) extracellular domain (19-740) was expressed CHO cells using a C-terminal Fc tag. The gene accession number is AF291820.

#### Alternative name(s)

ACEH.

#### **Formulation**

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 90mM glycine.

#### Storage and Stability

Store product at  $-70^{\circ}$ C. For optimal 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.

#### Scientific Background

The angiotensin converting enzyme 2 (ACE2) belongs to the angiotensin-converting enzyme family of dipeptidyl carboxydipeptidases (1). ACE2 cleaves the angiotensin I and angiotensin II to counterbalancing ACE activity in the regulation of renin angiotensin system (RAS) (1). ACE2 also has the role in regulation of cardiovascular, renal and reproductive systems (1). In addition, ACE2 is a functional receptor for the spike glycoprotein (S) of SARS-CoV, HCoV-NL63 (2), and SARS-CoV-2, which causes the pandemic of the respiratory diseases (COVID-19) around the world in 2020 (3).

#### References

- Imai Y, et al: Angiotensin-converting enzyme 2 (ACE2) in disease pathogenesis. Circ J. 2007, 74(3):405-10.
- Lan J, et al: Crystal structure of the 2019-nCov spike receptorbinding domain bound with the ACE2 receptor. bioRxiv. doi: https://doi.org/10.1101/2020.02.19.956235.
- Zhou P, et al: A pneumonia outbreak associated with a new coronavirus of probable bat origin. Nature. 2020, 579:270-89.

 Catalog #
 Aliquot Size

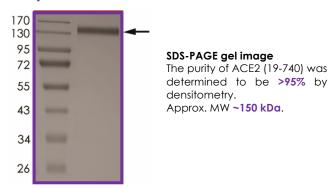
 A51C2-G341F-10
 10 μg

 A51C2-G341F-20
 20 μg

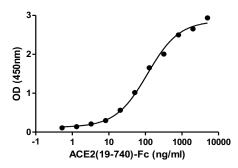
 A51C2-G341F-50
 50 μg

 A51C2-G341F-100
 100 μg

#### **Purity**



#### **Activity**



Measured by its binding ability in a functional ELISA. Immobilized 2019-nCoV spike protein RBD (His-tag) (C19SD-G241H) can bind to ACE2(19-740) protein (Fc tag).

# ACE2 (19-740) Protein

Recombinant human protein expressed in CHO cells

Catalog # A51C2-G341F Lot # O3814-4 Purity >95% Concentration  $0.5 \, \mu g/\mu l$ Stability 1yr at -70°C from date of shipment

Storage & Shipping

Store product at -70°C. For optimal storage and performance aliquot product into smaller quantities after centrifugation and store at recommended temperature. Avoid repeated handling and multiple freeze/thaw cycles.

Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: <a href="mailto:orders@signalchem.com">or IVD@signalchem.com</a> - www.signalchem.com

#### Revised date: 2022-09-06 Page 1 of 3

# SAFETY DATA SHEET

## **Article 1 - Product Identification**

#### Product Name: ACE2 (19-740) Protein

#### Catalog # A51C2-G341F

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. 110-13120 Vanier Place Street Address: Richmond, BC, V6V 2J2 City, Prov. Postal Code:

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

# **Article 2 - Hazard Identification**

- WHMIS Classification: Not WHMIS controlled.
- GHS classification: None.
- Hazard Pictograms: None.
- Signal words: None.
- Hazard statements: None.
- Precautionary statements: None.
- Other hazards: None known.

# Article 3 - Composition/Information on Ingredients

Chemical Characterization: Mixture.

**Description:** This product consists of the substances listed below.

Common name	Chemical name	CAS-No.	Concentration
Tris-HCl; Tris (hydroxymethyl) aminomethane	2 – Amino – 2 - (hydroxymethyl) propane -	1185-53-1	≤0.80%
hydrochloride	1, 3 - diol hydrochloride		
Glycine	Aminoethanoic acid	56-40-6	≤0.75%
Protein	-	No data available	≤0.05%

#### **Article 4 - First-aid Measures**

- General information: Consult a physician by providing the SDS.
- After inhalation: Breath in fresh air. If cannot breathe, give artificial respiration and consult a physician.
- After skin contact: Immediately wash with soap and plenty of water and rinse thoroughly. Consult a physician.
- 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: Not expected to present a significant ingestion hazard under anticipated conditions of normal use. If you feel unwell, seek medical advice.

#### 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.

# Revised date: 2022-09-06 Page 2 of 3

# SAFETY DATA SHEET

#### 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 safe 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 -70 °C. Keep container upright and tightly closed.

# **Article 8 - Exposure Controls/Personal Protection**

Components with limit monitoring values at workplace:

NA

- 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 available.
Boiling point/Boiling range: >100 °C.	Density: Not determined.
Flash point: > 100 °C.	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: Heat and moisture.
- Incompatible materials: Not determined.
- Hazardous decomposition products: Not determined.

# Revised date: 2022-09-06 Page 3 of 3

# SAFETY DATA SHEET

# **Article 11 - Toxicological Information**

- Acute toxicity: Not available.
- LD/LC50: Not available.
- Skin corrosion/irritation: Not available.
- Serious eye damage/eye irritation: Not available.
- 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: No data available
   Ingestion: No data available
   Skin: No data available
   Eyes: No data available
- Signs and Symptoms of Exposure: No data available
- Synergistic effects: Not available.

## Article 12 - Ecological Information

- Eco-toxicity: No data available.
- 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.