

DPP4 (29-766) Protein

Recombinant human protein expressed in CHO cells

Catalog # DP4CD-G441H

Lot # J3823-2

Product Description

Recombinant human DPP4 (29-766) Protein was expressed in CHO cells using a C-terminal His tag. The gene accession number is \underline{NM} 001935.

Alternative name(s)

ADABP; ADCP2; CD26; DPPIV; TP103

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.5, 300mM NaCl, 150mM imidazole.

Storage and Stability

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.

Scientific Background

The Middle East respiratory syndrome coronavirus (MERS-CoV) has caused the MERS outbreak that first originated in Saudi Arabia, but since has spread to other countries around the world (1). The spike glycoprotein (S) of coronavirus belongs to the type I transmembrane protein containing two subunits (2), S1 and S2, which is known to be the key component to bind with host cells through interaction with dipeptidyl peptidase 4 (DPP4) (3). MERS-CoV can recognize DPP4 through its receptor binding domain (RBD) of S1 to initiate viral infection in humans (2).

References

- Mohd H, et al: Middle East Respiratory Syndrome Coronavirus (MERS-CoV) origin and animal reservoir. Viral J. 2016, 13: 87.
- Lu G, et al: Molecular basis of binding between novel human coronavirus MERS-CoV and its receptor CD26. Nature. 2013, 500: 227-231.
- Al-Tawfiq et al: Middle East Respiratory Syndrome-Coronavirus (MERS-CoV) Infection. Emerging Infectious Diseases. 2014, 185-190.

 Catalog #
 Aliquot Size

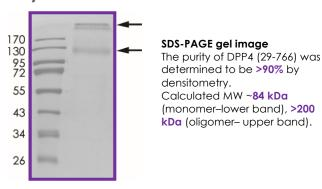
 DP4CD-G441H-10
 10 μg

 DP4CD-G441H-20
 20 μg

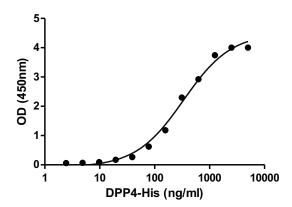
 DP4CD-G441H-50
 50 μg

 DP4CD-G441H-100
 100 μg

Purity



Activity



Binding ability measured in a functional ELISA. MERS Spike Protein RBD (MERSD-G241F) can bind to immobilized DPP4 (29-766) protein (DP4CD-G441H).

DPP4 (29-766) Protein

Recombinant human protein expressed in CHO cells

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: or IVD@signalchem.com - www.signalchem.com

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

SAFETY DATA SHEET

Article 1 - Product Identification

Product Name: DPP4 (29-766) Protein

Catalog # DP4CD-G441H

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
NaCl	Sodium chloride	7647-14-5	1.75%
Imidazole	1,3-Diaza-2,4-cyclopentadiene	288-32-4	≤1.02%
Sodium Phosphate, Dibasic	Sodium Phosphate, Dibasic	7782-85-6	1.34%
Protein	N/A	No data available	≤0.02%

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.