

p19INK4D Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # C33-30G

Lot # Q077-3

Product Description

Recombinant full-length human p19INK4D was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_001800](#).

Gene Aliases

CDKN2D, p19, INK4D, p19-INK4D

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°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.

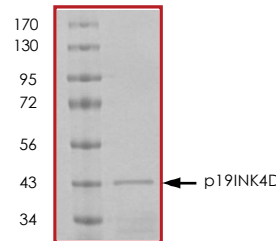
Scientific Background

p19INK4D protein is associated with CDK6 in several hematopoietic cell lines. It has been demonstrated that p19INK4D expression enhances cell survival under genotoxic conditions (1). By using p19INK4D-overexpressing cells, it has been demonstrated that p19INK4D expression correlates with the cellular resistance to UV treatment with increased DNA repair activity against UV-induced lesions. p19INK4D regulates protein network that integrate DNA repair, apoptosis and checkpoint mechanisms in order to maintain the genomic integrity (2).

References

1. Guan, K.L. et al: Isolation and characterization of p19INK4d, a p16-related inhibitor specific to CDK6 and CDK4. *Mol Biol Cell*. 1996; 7(1): 57-70.
2. Julieta, M. et al: Induction of p19INK4d in response to ultraviolet light improves DNA repair and confers resistance to apoptosis in neuroblastoma cells. *Oncogene*, 2005; 24: 4065-4080

Purity



The purity of p19INK4D was determined to be **>90%** by densitometry. Approx. MW **44kDa**.

p19INK4D Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number C33-30G

Specific Lot Number Q077-3

Purity >90%

Concentration 0.2 µg/µl

Stability 1 yr at -70°C from date of shipment

Storage & Shipping 1 yr at -70°C from date of shipment

Store product at -70°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. 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: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

MATERIAL SAFETY DATA SHEET

Article 1 - Product Identification and Use

Product Name: p19INK4D Protein

Catalog # C33-30G

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 Pharmaceuticals Inc.
Street Address: 550-5600 Parkwood Way
City, Prov. Postal Code: Richmond, BC, V6V 2M2
Fax: 604-232-4601
EMERGENCY PHONE: 604-232-4600

Article 2 - Hazardous Ingredients

NOT AVAILABLE. We are not aware of any hazards associated with this product or its ingredients, but the chemical, physical, and toxicological properties of this product have not been investigated thoroughly. Observe normal laboratory precautions.

Article 3 - Physical Data

This product consists of purified protein in Tris-HCl buffer shipped on dry ice. The physical properties of this product have not been investigated thoroughly.

Article 4 - Fire and Explosion Hazard

NOT APPLICABLE

Article 5 - Reactivity Data

NOT APPLICABLE

Article 6 - Toxicologically Data

May be harmful by inhalation, ingestion, or skin absorption. The toxicological properties of this product have not been investigated thoroughly. Exercise due caution.

Article 7 - Preventative Measures

Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact with eyes, skin or clothing.

*****MULTIPLE COMPONENT SPILL OR LEAK PROCEDURES*****

- Wear protective equipment.
 - Absorb on sand or vermiculite and place in closed containers for disposal.
 - Observe all federal, state and local environmental regulations.
-

Article 8 - First Aid Measures

- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
 - In case of skin contact, flush with copious amounts of water for at least 15 minutes. Remove contaminated clothing and shoes. If a rash or other irritation develops, call a physician.
 - If inhaled, remove to fresh air. If breathing becomes difficult, call a physician.
 - In case of eye contact, flush with copious amounts of water for at least 15 minutes while separating the eyelids with fingers. Call a physician.
-

Article 9 - Preparation

Prepared by: Jun Yan

Phone#: 1-866-954-6273

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 Catalogue for additional terms and conditions of sale.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.