

Datasheet

V. 2.3

Cytoskeleton, Inc.

Ultra-Precise HRV 3C Protease (human rhinovirus type 14 3C Protease) (recombinant-GST tagged) Cat. # CS-UPP01 Lot # 013 10 x 100 Units Upon arrival store at 4°C (desiccated) See datasheet for storage after reconstitution

Background Information

The Ultra-Precise HRV 3C Protease (UP protease) is a fusion protein with a glutathione S-transferase affinity tag (GST) and human rhinovirus (HRV) type 14 3C protease commonly used in molecular biology and protein purification techniques. The UP protease specifically cleaves the specific amino acid sequence Leu-Glu-Val-Leu-Phe-Gln/Gly-Pro (LEVLFQ/GP), cleaving between the GIn and Gly residues. UP protease can maintain robust cleavage activity at 4°C, allowing for improved target protein stability. The GST-tag of UP protease allows for the simultaneous cleavage of a targeted GST-fusion protein substrate and the protease removal when cleaving on Glutathione Sepharose Resin. The cleaved protein substrate can be collected in the flow-through while leaving GST and UP protease bound to the resin. The UP protease is identical in sequence and activity to the popular PreScission protease. The UP protease cleavage site can be found in fusion proteins produced from pGEX-6P vectors.

Material

The recombinant UP protease protein was produced in a bacterial expression system. The recombinant protein has an N-terminal GST-tag and has an approximate molecular weight of 46 kDa. One unit is the amount of enzyme capable of cleaving >90% of 100 μ g of fusion protein in 16 hours at 4°C. Protease specific activity is 800-1100 units/mg. The protein is supplied as a white solid.

Storage and Reconstitution

The recommended storage conditions for the lyophilized material are desiccated at 4°C. Under these conditions, the protein is stable for one year. Lyophilized protein can also be stored desiccated at -70° C.

Before reconstitution, briefly centrifuge to collect the product at the bottom of the tube. The protein should be reconstituted by adding 100 μ l of cleavage buffer for 1 U/ μ l. When reconstituted, the protein will be in the following buffer: 50 mM Tris pH 7.0, 150 mM NaCl, 1 mM EDTA, 1 mM DTT, 1.5% sucrose, and 0.2% Dextran. Stable at 4°C for 1-2 weeks.

Figure 1. Purity analysis of UP Protease residues

Legend: A 10 µg sample of UP protease purity is

determined by scanning densitometry of Coomassie

Blue-stained protein on a 4-20% polyacrylamide

gradient gel. The UP protein was determined to be

90% pure. (see Figure 1). Protein quantitation was

performed using the Precision Red Protein Assay

Reagent (Cat. # ADV02). Molecular weight markers

Cleavage Buffer

The following cleavage buffer needs to be prepared before working with UP Protease.

50 mM Tris-HCl pH 7.0 150 mM NaCl 1 mM EDTA 1 mM DTT

Note: Digestion may be improved by adding 0.01% detergents (Triton X-100, Tween-20, and NP-40). Concentrations up to 1% of detergents do not inhibit protease activities. The presence of Zrf⁺ has been found to reduce UP protease activity.

Small-scale cleavage optimization of GST-Fusion proteins.

The cleavage activity of UP protease should be optimized on a small scale before scaling up. The amount of UP protease and digestion length for the fusion protein of interest may vary.

Reagents

- 1. Ultra-Precise HRV 3C Protease (Cat. # CS-UPP01)
- 100 µg of GST-Fusion Protein
- 3. Cleavage Buffer at 4°C

Method

- 1. Resuspend 100U of UP Protease with 100 μl of cleavage buffer.
- Ensure the GST-Fusion protein has been dialyzed or resuspended in cleavage buffer.
- 3. Add 100 μg to each of the 4 reaction tubes labeled 2U, 1U, 0.5U, and 0U respectively.
- Add 2 μl (2 U), 1 μl (1U), 0.5 μl (0.5U), and no UP protease as a control to the corresponding reaction tubes. Higher or lower dilution reactions can be tested if desired.
- Incubate for 16hrs at 4°C.
- After 16hrs, add 20 μl of reaction to 20 μl of 2X SDS-PAGE loading buffer.
- Determine cleavage activity by running samples on an SDS -PAGE gel.
- Reactions can be scaled up after determining ideal digestion conditions.
- 9. The time for digestion can also be determined with this method with a fixed unit of UP Protease and removing 20 μ I samples at different time points.

cytoskeleton.com

are from Invitrogen (Mark 12).

Phone: (303) 322.2254 Fax: (303) 322.2257 Customer Service: cserve@cytoskeleton.com Technical Support: tservice@cytoskeleton.com



Figure 2. Cleavage of fusion-protein with UP protease Legend: 100 μ g of GST- α TAT1 protein was cleaved with UP protease according to the "Small scale cleavage test" conditions in this datasheet. A 20 ug sample of each condition was run on SDS-PAGE, and cleavage efficiency was determined by densitometric analysis of Coomassie blue stained protein bands. Different units of UP were used per reaction (2 units, 1 unit, and 0.5 unit) to determine the amount of enzyme needed for digestion. Molecular weight markers are from Invitrogen (Mark 12).

On-column Cleavage Method

On-resin cleavage of the GST-fusion protein of interest with UP protease is recommended, as the cleaved protein of interest can be separated from GST and UP protease.

- The recombinant GST-fusion protein of interest can be bound to glutathione resin from cell lysate after the chosen cell lysis method.
- The glutathione resin can be washed with 5-10 bed volumes of chosen wash buffer to remove protein impurities.
- 3. Exchange glutathione resin with cleavage buffer and allow the resin to settle. Ensure at least one equal bed volume of cleavage buffer rests above the resin bed. Add 1 to 2U of UP protease per 100 μg of GST-fusion protein and incubate the reaction for 16hrs at 4°C. The reaction can be rotated/ shaken slowly overnight to help ensure complete digestion, but specific proteins may precipitate with rotation/shaking versus allowing the reaction to sit overnight with no agitation.
- 4. The protein of interest can be eluted and collected.
- 5. The UP protease will remain bound to the glutathione resin.