

Product Information

Thrombin CleanCleave™ Kit

Catalog Number **RECOMT**
 Storage Temperature 2–8 °C

TECHNICAL BULLETIN

Product Description

The Thrombin CleanCleave™ Kit contains a 50% (v/v) suspension of thrombin-agarose produced by immobilizing bovine thrombin (molecular mass ~37 kDa) and is designed for cleavage of recombinant fusion proteins. This format circumvents the need for removal of thrombin after cleavage by chromatographic techniques. The ligand density used allows for fast and efficient cleavage; the resin can be reused multiple times with only minimal loss in cleavage efficiency.

Thrombin-agarose can be used to cleave any fusion protein containing a thrombin cleavage site. The optimal cleavage sites for thrombin are:¹

1. P4-P3-Pro-Arg/Lys • P1'-P2'
 P4 and P3 are hydrophobic residues, P1' and P2' are non-acidic residues, and Arg/Lys • P1' is the scissile bond.
2. P2-Arg/Lys • P1'
 Where P2 or P1' is glycine and Arg/Lys • P1' is the scissile bond.

Cleavage of the fusion protein is performed in a batch format and removal of the resin is accomplished either by centrifugation or by collection of the resin in a column. The cleavage time may range from 2 hours to overnight depending upon reaction conditions, temperature, amount of resin used, and fusion protein.

Thrombin-agarose has a wide range of activity.² The optimal pH range for cleavage of a fusion protein is between 8–9. Thrombin-agarose is active at a pH as low as 5.0; however, cleavage rates are reduced by more than 40% when the pH of the reaction is dropped below 7, see Figure 1.

Thrombin-agarose can be used over a variety of temperatures from 4–37 °C. Cleavage rates at 4 °C are significantly reduced (80%) compare to rates at room temperature. A minimal increase in cleavage rate is observed at 37 °C compared to 25 °C.

The resin can also be used over a wide ionic strength range and in the presence of number of buffer additives allowing for compatibility with an assortment of recombinant protein expression strategies. Cleavage was significant from 0–500 mM NaCl, with only a 20% activity change in cleavage over this NaCl range. Thrombin-agarose has been found to function well in the presence of 0.1% TRITON® X-100, 1 M urea, and 5 mM EDTA. This is by no means a comprehensive list.

Thrombin contains disulfide bonds, so reducing agents such as dithiothreitol and 2-mercaptoethanol should be avoided.

Resin Characteristics

Matrix: 4% Beaded agarose

Ligand immobilized: ≥0.15 mg/ml resin

Molecular weight exclusion limit: 2×10^7

Particle size: 45–165 µm

Maximum linear flow rate: ~11.5 cm/hr

Cleavage rate of synthetic substrate: 10 µmoles of *N*-benzoyl-Phe-Val-Arg-*p*-nitroanilide (Catalog Number B7632) per µmole of bound thrombin per second.

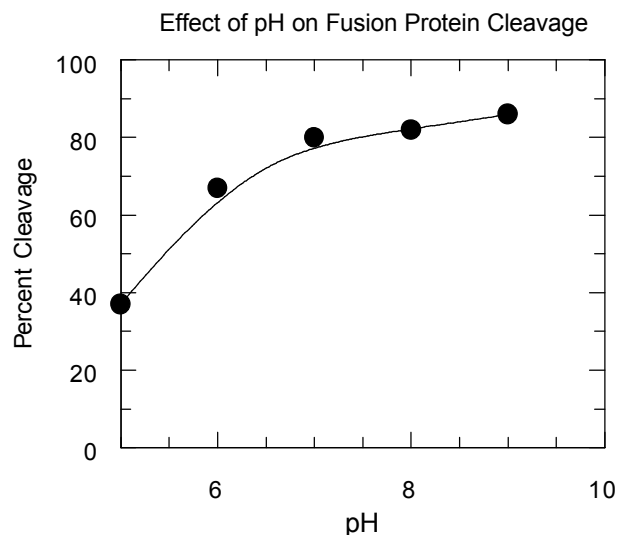
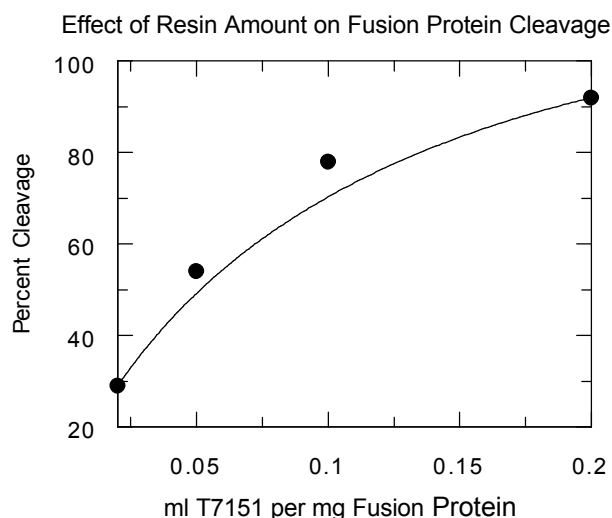
Cleavage rate of fusion protein: 200 µl of a 50% suspension of resin will cleave >85% of 1 mg of fusion protein in 1 ml in 4 hours at room temperature, see Figure 2.

Stability of linker: No detectable leakage of thrombin from thrombin-agarose after 1 month at 2–8 °C.

Detection limit: <5 picomoles/ml packed gel

Components

Thrombin-Agarose	1 ml
50% suspension in 50% glycerol, 20 mM Tris-HCl, pH 8.2. Furnished as ml packed gel. Catalog Number T7151	
10× Cleavage Buffer	10 ml
500 mM Tris-HCl, pH 8.0, 100 mM CaCl ₂ Catalog Number T9685	
Disposable column	1 each
Catalog Number C4969	

Figure 1.**Figure 2.****Precautions and Disclaimer**

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation InstructionsSample Preparation

Samples of fusion proteins should be purified prior to cleavage. Samples having a high viscosity due to the presence of chromosomal DNA or RNA will clog the resin. Purification is best accomplished by passing the sample over an affinity column specific for the fusion protein. The fusion protein should be exchanged into 50 mM Tris-HCl, pH 8.0, with 10 mM CaCl₂. Samples that contain particulates should either be centrifuged or filtered. Crude mixtures may reduce the reusability and reproducibility of the thrombin agarose. Samples containing proteases should be used with caution as these samples may also reduce the performance of the resin. Protease inhibitors may reduce the performance of the resin and should be avoided.

Storage/Stability

Store the kit at 2–8 °C.

ProcedureFusion Protein Cleavage Procedure

These guidelines are for cleaving of 1 mg of fusion protein in a final volume of 1 ml. Determining optimal cleavage conditions, temperature, the length of incubation, and the stability of the fusion protein, should be done on pilot experiments to achieve maximum recovery prior to large scale reactions.

Washing Resin

1. Thoroughly resuspend the thrombin-agarose resin to make a homogeneous slurry.
2. Remove a 100 µl aliquot of a 50% (v/v) suspension of resin and gently spin in a microcentrifuge at 500 × g (2,500 rpm) to pellet the resin. Remove the supernatant.
3. Add 500 µl of 1× Cleavage Buffer and gently resuspend by inverting the tube. Centrifuge at 500 × g and remove the supernatant.
4. Repeat step 3.

Cleavage Reactions

1. Add 100 µl of 10× Cleavage Buffer to the centrifuged beads and gently resuspend the beads.
2. Add 1 mg of fusion protein and bring to a final volume of ~1 ml with water.
3. Incubate the cleavage reaction at room temperature or other appropriate temperature with gentle agitation to keep beads suspended.
4. Remove aliquots at 1, 2, 4, 6, and 24 hours, and gently centrifuge to remove the resin and analyze supernatant for cleavage. Analysis of cleavage reaction may be done by SDS-PAGE.

Recovery of Fusion Protein

Select one of the two procedures for elution.

A. Centrifugation Recovery

1. Centrifuge the mixture for 5 minutes at $500 \times g$. Remove the supernatant containing the cleaved fusion protein.
2. To increase recovery, resuspend the resin with an equal volume of $1\times$ Cleavage Buffer, centrifuge the mixture, remove supernatant, and add to the first supernatant.
Note: The resin may be centrifuged briefly at high g forces ($10,000 \times g$) with minimal loss in cleavage efficiency. In order to prolong the effectiveness of the resin; however, a 5 minute centrifugation at $500 \times g$ is recommended.

B. Column Recovery

1. Transfer the cleavage reaction solution into the column.
2. Allow the resin to settle. Drain and collect the eluent.
3. Rinse the resin with 5 bed volumes of $1\times$ Cleavage Buffer and collect the eluent.
4. Combine eluents.

Resin Regeneration

1. Rinse the resin with an equal volume of 50 mM Tris-HCl, pH 8.0, with 500 mM NaCl.
2. Rinse the resin with an equal volume of 50 mM Tris-HCl, pH 8.0.
3. Repeat step 2 twice.
4. Store as a 50% suspension in 50% glycerol, 20 mM Tris-HCl, pH 8.0, at $2-8^\circ\text{C}$.
Note: Thrombin should be stored in the absence of NaCl or other salts to prolong the life of the resin.

The resin can be used a minimum of five times with little loss of effectiveness. Reusability will depend upon cleavage conditions, incubation time, temperature, sample preparation, the use of chaotropic or other extreme conditions, and proper storage condition.

Results

Troubleshooting

1. No cleavage seen:
 - Check sequence to verify there is an appropriate thrombin binding site.
 - Use optimal cleavage conditions.
2. Incomplete cleavage:
 - Increase the resin to fusion protein ratio and increase the incubation time to overnight.
 - Use optimal pH, salt concentration, and increase the temperature to 25°C or higher if necessary.
 - The addition of 1 M urea or other mild denaturants may be useful.
3. Secondary cleavage:
 - A second thrombin site may exist.
 - Reduce the incubation time and the resin to fusion protein ratio to minimize secondary cleavage.
 - Use suboptimal cleavage conditions to slow the cleavage rate down.
 - Reduce the temperature of the cleavage reaction to $2-8^\circ\text{C}$.

References

1. Chang, J.-Y., Eur. J. Biochem., **151**, 217 (1985).
2. Di Cera, E. *et al.*, Biochemistry, **30**, 7913-7924 (1991).

CleanCleave is a trademark of Sigma-Aldrich® Biotechnology LP and Sigma-Aldrich Co.
TRITON is a registered trademark of Dow Chemical Co.

MAM 07/10-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.