

Product Information

CellLytic[™] B 2X

Product Code **B 7310**
Store at Room Temperature

TECHNICAL BULLETIN

Product Description

CellLytic[™] B 2X is used for the lysis of bacterial cells for the purification of recombinant and wild type proteins. CellLytic B 2X contains 40 mM Tris-HCl, pH 8.0 and a proprietary, non-denaturing formulation of zwitterionic detergents. CellLytic B 2X is different from CellLytic B in that it provides a more concentrated protein solution. There is no need for special equipment to disrupt cells such as a sonicator or French press. This fast one-step method obtains higher yields than sonication, lysozyme treatment, or other commercially available lysis solutions. CellLytic B 2X can be used to extract soluble proteins and can also be used to wash away cell debris from inclusion bodies to yield nearly pure protein. CellLytic B 2X does not solubilize inclusion bodies.

CellLytic B 2X is optimized for the lysis of *E. coli* strain BL21. However, it also works well for DH5 α [™], JM109, and other similar bacterial cells. Intact fusion proteins have been successfully purified from CellLytic B lysates of BL21 *E. coli* cells expressing histidine-tagged and FLAG-tagged proteins using HIS-Select[™] and Anti-FLAG[®] M2 purification resins, respectively. The CellLytic B 2X reagent is also compatible with affinity purification of other fusion proteins.

Reagents and Equipment Required but Not Provided

(Product Codes have been given where appropriate)

- Lysozyme Solution (Product Code L 3790)
- Benzonase[®] (Product Code E 1014)
- Deionized or molecular biology grade water (Product Code W 4502), for the dilution of the CellLytic B reagent
- HIS-Select Nickel Affinity Gel (Product Code P 6611) or other fusion protein purification system
- Protease Inhibitor Cocktails for:
 - Bacterial Cells (Product Code P 8465)
 - Histidine-tagged Proteins (Product Code P 8849)
- CellLytic IB (Product Code C 5236)
- Appropriate centrifuge tubes
- Centrifuge

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.

It is recommended that the entire technical bulletin be read prior to use, especially the reagent compatibility chart.

Preparation Instructions

CellLytic B 2X is supplied as a ready-to-use solution. Additional reagents may be added to suit particular needs. These include salts, protease inhibitors, EDTA, and reducing agents such as dithiothreitol or 2-mercaptoethanol.

Storage/Stability

Store the product at room temperature. CellLytic B 2X, as supplied, is stable for at least 2 years at room temperature.

Procedures

A. Trial Scale Extraction

A small-scale trial extraction should be performed to determine the fraction in which the protein of interest will be found. If this has already been determined, proceed to the Large Scale Extraction section.

1. A five ml culture of the bacterial strain containing the recombinant protein should be grown under the appropriate conditions for expression.
2. Use 1.5 ml of the bacterial culture with an OD₆₀₀ of 0.5-1.0 and centrifuge the cells at full speed for 2 minutes.
4. Remove the spent medium and resuspend the cell pellet in 0.2 ml of CellLytic B 2X.
5. Briefly vortex the solution to resuspend the cell pellet and mix for 5-10 minutes to ensure full extraction of the soluble proteins.

6. Centrifuge the cell lysate at full speed for 5 minutes to pellet any insoluble material.
7. Carefully remove the soluble protein fraction from the cell debris. Additional extractions may be performed if required; however, this will result in a more dilute soluble protein sample.
8. Analyze the supernatant and the insoluble fraction by SDS-PAGE and/or Western blot to determine which fraction contains the protein of interest. For SDS-PAGE, it is recommended that 5-15 μ l of each sample be applied to the gel.

Note: If the protein of interest is not found in the soluble portion, it has likely formed inclusion bodies. For the purification/solubilization of inclusion bodies see the Inclusion Body Purification Procedure, Section C.

B. Large Scale Extraction

This procedure is designed for 1 gram of wet cell paste. This is roughly equivalent to a 250 ml bacterial culture with an OD₆₀₀ of ~2.0. In order to extract the maximum amount of soluble protein, the CelLytic B 2X to cell mass ratio should be 5 ml per gram of wet cell paste. Using less CelLytic B 2X will give a more concentrated solution, but a smaller amount of total protein will be extracted. Using more CelLytic B 2X will not extract any more protein; it will only serve to provide a more dilute protein solution.

1. Collect the bacterial cells that express the protein of interest by centrifuging at 5,000 x *g* for 10 minutes.
2. Carefully remove the spent medium from the cell pellet. The cell pellet may be frozen or used fresh. A frozen cell pellet will give a slightly higher yield of protein.
3. Add CelLytic B 2X at a ratio of 5 ml per gram of cell paste. Mix well to completely resuspend the cells. The following optional reagents may also be added:
 - a. Lysozyme (final concentration of 0.2 mg/ml) to enhance cell lysis.
 - b. Benzonase (final amount of 50 units/ml) to decrease the viscosity of the solution.
 - c. Protease inhibitors to prevent proteolytic degradation.
4. Incubate the extraction suspension with shaking at room temperature for 10-15 minutes to fully extract the soluble proteins from the cells.

5. After the cells have been extracted, centrifuge the extract at 16,000 x *g* for 10 minutes to pellet the insoluble material.
6. Carefully remove the supernatant containing the soluble protein fraction. Another round of extraction will yield more soluble protein if required; however, this will result in a more dilute soluble protein sample.
7. Analyze the supernatant and insoluble fraction by SDS-PAGE and/or Western blot to determine which fraction contains the protein of interest. For SDS-PAGE, it is recommended that 5-15 μ l of each sample be applied to the gel.

Note: CelLytic B will **not** solubilize inclusion bodies. For purification of inclusion bodies, see the Inclusion Body Purification Procedure, Section C.

C. Inclusion Body Purification

1. Resuspend the cell pellet from the first extraction (Section A, step 7 or Section B, step 6) in a volume of CelLytic B 2X solution equal to that was used for the extraction. Vortex for 1-2 minutes to completely resuspend the cell debris. Add Lysozyme Solution (Product Code L 3790) to a final concentration of 0.2 mg/ml.
2. Incubate at room temperature for 5-10 minutes to allow the lysozyme to fragment the cell wall.
3. Dilute CelLytic B 2X 20-fold by mixing 95 ml of deionized water for every 5 ml of CelLytic B.
4. Add 30 ml of the 20-fold diluted CelLytic B 2X per gram of paste from the original extraction to the tube and mix thoroughly.
5. Centrifuge at full speed for 5 minutes to pellet the cell debris again. Save the supernatant for analysis.
6. Resuspend the pellet in a volume of 20-fold diluted CelLytic B 2X equal to that was used in step 4 and vortex to completely resuspend any remaining insoluble material.
7. Centrifuge at full speed for 5 minutes to pellet the cell debris. Save the supernatant for analysis. Steps 5 and 6 may be repeated a number of times to completely remove any remaining soluble proteins and cell wall material from the inclusion bodies. This wash step should be optimized for the specific protein of interest.

8. Resuspend the washed inclusion bodies (pellet from step 6) in an equal volume of deionized water or a buffer of choice.
9. Analyze all of the saved supernatants and the insoluble fraction by SDS-PAGE and/or Western blot. For SDS-PAGE, it is recommended that 5-15 μ l of each sample be applied to the gel.

Note: Alternatively, inclusion bodies can be solubilized in CellLytic IB (Product Code C 5236).

Reagent Compatibility Chart

Reagent	Effect	Comments
Chelating agents (EDTA, EGTA)	Strips metal ions from IMAC resins and chelates essential Mg^{2+}	EDTA is not compatible with the HIS-Select line of products. It will chelate metal ions from the affinity gel. Also, addition of EDTA to the original cell lysis mixture will chelate metal ions essential for endonuclease activity, which will result in a thick, viscous solution.
Protease Inhibitors	Prevent protein degradation	Protease inhibitors may be added to the bacterial cell culture extraction, if desired.
2-mercaptoethanol and dithiothreitol	Reducing Agents	Can be used at low levels for downstream application to HIS-Select products; should not be used for FLAG or glutathione resins.

Troubleshooting Guide

Problem	Cause	Solution
Lower than expected protein levels	Cells are not completely lysed.	Freeze/thaw cells to increase cellular breakage. Addition of lysozyme (final concentration of 0.2 mg/ml) will aid in protein extraction.
	Sample viscosity is too high.	Addition of Benzonase (final activity of 50 units/ml) will reduce sample viscosity and aid in recovery of soluble extract.
	Target protein degraded.	Addition of protease inhibitors may help reduce target protein degradation. See Reagents and Equipment Required but Not Provided for recommended protease inhibitor cocktails.
	Expression level may be too low.	<ul style="list-style-type: none"> Add more inducing agent. Induce for a longer time period. Check the construct. Use another bacterial cell line.
	Protein of interest may be insoluble.	Check pellet to ensure protein of interest has not formed inclusion bodies.
	Not enough CellLytic B added to cells for full extraction.	Use more of the reagent per amount of cells. However, it is not beneficial to add more than 10 ml per gram of wet cell paste

Benzonase is a registered trademark of Merck KGaA.
DH5 α is a trademark of Invitrogen.

JP/MAM 03/05-1

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.