

Dr. Gold's Bacterial Protein Extraction Protocol For the Extraction of Expressed Proteins

Introduction

Dr. Gold's Bacterial Protein Extraction Reagent is a powerful, next-generation solution designed for efficient extraction of expressed proteins from bacterial cells—now reengineered for improved performance and cost-effectiveness. Built on the foundation of our trusted Bacterial Cell Lysis Buffer, this ready-to-use formulation offers enhanced yield and reproducibility without the need for harsh mechanical disruption.

The reagent includes lysozyme, nucleases, and a specially optimized salt solution to promote full membrane disruption and maximize the recovery of soluble proteins.

Note: The Salt Solution is not strictly required for lysis, but our research has shown they significantly enhance protein solubility and yield.

Materials

- Dr. Gold's Bacterial Protein Extraction Reagent
- GB Enzyme Mix
- Salt Solution
- Pipet
- Centrifuge tubes
- Microcentrifuge tubes
- Centrifuge
- Shaker incubator
- Incubator (22°C)
- UV/Vis Spectrophotometer

Storage and Handling

- This product may be shipped with blue ice. Dr. Gold's Bacterial Protein Extraction Reagent and Salt Solution should be stored at 4°C, GB Enzyme Mix should be stored at -20°C immediately upon arrival. When stored under the recommended conditions and handled correctly, these products should be stable for at least 1 year from the date of receipt.
- Thaw the **GB Enzyme Mixture** on ice. Gently flick the tube to resuspend any settled material.

Method

Note: This protocol assumes you have a culture of *E. coli* cells expressing your protein of interest. The culture should be properly induced and grown under optimized conditions. For example, *E. coli* cells carrying a plasmid with an IPTG-inducible promoter (e.g., lac) may be induced with 0.1M IPTG and grown overnight at 30°C.

1. **Measure cell density**

Measure the optical density at 600 nm (A_{600}) of the *E. coli* culture. You will need this to determine the amount of required lysis buffer on step 3.

2. **Harvest the cells**

Centrifuge the culture at **3,500 × g for 15 minutes** (or 10,000 × g for 5 minutes) to pellet the cells. *Note:* 15 minutes at 3,500 × g is generally ideal if your centrifuge rotor has a 12 cm path length.

3. **Resuspend in lysis buffer**

Discard the supernatant and gently resuspend the pellet in **pre-warmed (22°C) lysis buffer**. If desired, you may add protease inhibitors to the lysis buffer prior to resuspending.

Note: For Protease inhibitors, we recommend using ProBlock™ Gold Bacterial Protease Inhibitor Cocktail (GoldBio Catalog # GB-330)

Use **1.3 mL of lysis buffer per 10 mL of culture at $A_{600} = 2.0$** . Adjust proportionally for different cell densities and volumes.

Important: Do NOT vortex. Resuspend by gentle pipetting or swirling.

4. **Add enzyme mixture**

Thaw the **GB Enzyme Mixture** on ice. Gently flick the tube to resuspend any settled material.

Add **1/100 volume** of enzyme mix to the resuspended cells. Mix gently.

5. **Incubate for enzymatic lysis**

Incubate at **22°C for 30 minutes**, gently agitating every 10 minutes.

Observation: Some debris may settle as the enzymes digest cell walls, DNA, and RNA.

6. Add salt solution

Add **1/10 volume** of the included **Salt Solution** to the lysate.

The solution should become more translucent, though some opaqueness may remain. Place the mixture on **ice** for 5-10 minutes.

Note: If analyzing lysate via SDS-Page, remove a 0.5 mL aliquot into a microcentrifuge tube and process separately for soluble/insoluble fraction analysis (See instructions for Optional Analysis via SDS-Page below).

7. Clarify lysate

Centrifuge the mixture at **10,000 × g for 10 minutes at 4°C** to pellet cell debris.

8. Harvest supernatant

Carefully transfer the **clarified supernatant** (soluble protein fraction) for downstream analysis or purification.

Analysis via SDS-PAGE (Optional):

1. For the **lysate (supernatant)**: After centrifugation according to **Step 8**, remove supernatant and mix directly with SDS-Page loading buffer for evaluation.
2. For the **pellet (insoluble fraction)**: Resuspend pellet in **50 µl of SDS-Page loading buffer**, heat at **100°C for 3 minutes**, then centrifuge at **10,000 × g for 5 minutes**.
3. Analyze both supernatants by SDS-PAGE to evaluate soluble vs. insoluble protein.

Associated Products

- Dr. Gold's Bacterial Protein Extraction Reagent (GoldBio Catalog # GB-500)
- ProBlock™ Gold Bacterial Protease Inhibitor Cocktail (GoldBio Catalog # GB-330)