

In-gel Digestion of Proteins

(Modified from University of Wisconsin – Biotechnology Center protocol)

Gel fragment preparation

- Excise protein bands. Cut each into 1 mm pieces. Place into a low-binding centrifuge tube.
- Wash gel pieces twice with copious dH₂O (~200 µl) for 30 seconds each to wash out acetic acid.

Destaining

- **Coomassie Blue** – ([Coomassie Brilliant Blue G-250](#) or [Coomassie Brilliant Blue, R-250](#))
Destain two times for 5 minutes or until colorless with 200 µl 100mM (NH₄)HCO₃/50% Methanol (discard supernatants). Dehydrate for 5 minutes with 200 µl 25mM (NH₄)HCO₃/50% Acetonitrile, then once more for 30 seconds in 100% Acetonitrile. Remove the solutions and discard. The gel pieces shrink and become white.
- **Non-Destructive Silver staining** – Destain twice for 2 minutes or until colorless with 200 µl of freshly prepared 1:1 solution of 100mM Sodium Thiosulfate [Na₂S₂O₃] and 30mM Potassium Ferricyanide [K₃Fe(CN)₆]. Stop the reaction and wash out silver ions twice for 2 minutes with 500 µl of Millipore water. Dehydrate for 5 minutes with 200 µl 25mM (NH₄)HCO₃/50% Acetonitrile, then once more for 30 seconds in 100% Acetonitrile. The gel pieces shrink and become white.
- **Sypro Ruby staining** – No destaining necessary. Dehydrate for 5 minutes with 200 µl 25mM (NH₄)HCO₃/50% Acetonitrile, then once more for 30 seconds in 100% Acetonitrile. Remove the solutions and discard. The gel pieces shrink and become white.
 - Dry gel particles for 2 minutes in a vacuum centrifuge.

Reduction and Alkylation

1. Prepare fresh:
 - 25mM Dithiothreitol ([DTT; GoldBio Catalog # DTT](#)) in 25mM Ammonium Bicarbonate ((NH₄)HCO₃), and
 - 55mM Iodoacetamide in 25mM Ammonium Bicarbonate.
2. Rehydrate gel pieces in 50 µl of prepared 25mM DTT. Reduce the proteins for 20 minutes at 56°C.
3. Cool the gel fragments to room temperature, pipet off any residual liquid and add 50 µl of 55mM Iodoacetamide. Alkylate the proteins for 20-30 minutes at room temperature in the dark.
4. Remove liquid and wash gel pieces with ≥20 volumes of dH₂O water (~400 µl) for 30 seconds to remove residual Iodoacetamide. Dehydrate gel pieces for 5 minutes with 200 µl 25mM (NH₄)HCO₃/50% Acetonitrile.

Gold Biotechnology

St. Louis, MO

Web: www.goldbio.com

Ph: (314) 890-8778

email: contactgoldbio86@goldbio.com

5. Remove liquid and dehydrate gel pieces again for 30 seconds in 100% Acetonitrile.
6. Dry gel particles for 2 minutes in a vacuum centrifuge.

Digestion

1. Prepare fresh Trypsin solution (10 ng/ μ L in 25mM Ammonium Bicarbonate/3% Acetonitrile, pH ~8.5).
2. Rehydrate with 40 μ L Trypsin solution and incubate for 4-5 hours (or overnight, if desired) at 37°C.
3. Overlay the rehydrated gel particles (if necessary) with a minimum amount of 25mM (NH₄)HCO₃ to keep them immersed throughout the digestion.

Peptide Recovery

1. Extract digested peptides with 50 μ L dH₂O/1% TFA (Trifluoroacetic Acid) by vortexing for 10 minutes at room temperature. Transfer solution to a new low-binding tube.
2. Perform an additional extraction with 80 μ L of 70% acetonitrile/25% H₂O/5% TFA.
3. Dry peptide solution completely in a vacuum centrifuge (1-2 hours).
4. Reconstitute peptides in 20 μ L dH₂O/0.1% TFA by incubating for 5 minutes at room temperature with intermittent vortexing.
5. Sample is ready for LC-based analysis.

References

Gharahdaghi, F., Weinberg, C. R., Meagher, D. A., Imai, B. S., & Mische, S. M. (1999). Mass spectrometric identification of proteins from silver-stained polyacrylamide gel: a method for the removal of silver ions to enhance sensitivity. *Electrophoresis*, 20, 601-605.

Jiménez, C. R., Huang, L., Qiu, Y., & Burlingame, A. L. (1998). In-Gel Digestion of Proteins for MALDI-MS Fingerprint Mapping. *Current Protocols in Protein Science*, 16-4.

Katayama, H., Nagasu, T., & Oda, Y. (2001). Improvement of in-gel digestion protocol for peptide mass fingerprinting by matrix-assisted laser desorption/ionization time-of-flight mass spectrometry. *Rapid Communications in Mass Spectrometry*, 15(16), 1416-1421.