

HaloTag[®] Mammalian Protein Detection and Purification Systems

Quick Protocol

Instructions for Use of Products **G6790 and G6795**

HaloTag[®] Purification—Quick Purification Protocol from 2×10^8 cells

Lyse

1. Resuspend the cell pellet in 5ml of HaloTag[®] Purification Buffer
2. Add 100 μ l of 50X Protease Inhibitor Cocktail.
3. Sonicate on ice (avoid overheating as this will inhibit binding).
Note: For other lysis methods, refer to Technical Manual TM348.
4. Harvest cell lysate at 4°C (10,000 \times g for 15 minutes); collect supernatant.

Equilibrate Resin

5. Transfer 600 μ l of HaloLink[™] Resin slurry to a tube.
6. Centrifuge at 1,500 \times g for 5 minutes; discard the supernatant.
7. Wash the resin five times:
 - a. Add 5ml of HaloTag[®] Purification Buffer; mix for 5 minutes.
 - b. Centrifuge at 1,500 \times g for 5 minutes; discard the supernatant.

Bind

8. Add the cell lysate to the equilibrated resin.
9. Incubate for 90 minutes at room temperature (22–25°C) with constant mixing.
10. Centrifuge at 1,500 \times g for 5 minutes; remove supernatant, save as sample flowthrough.

Wash

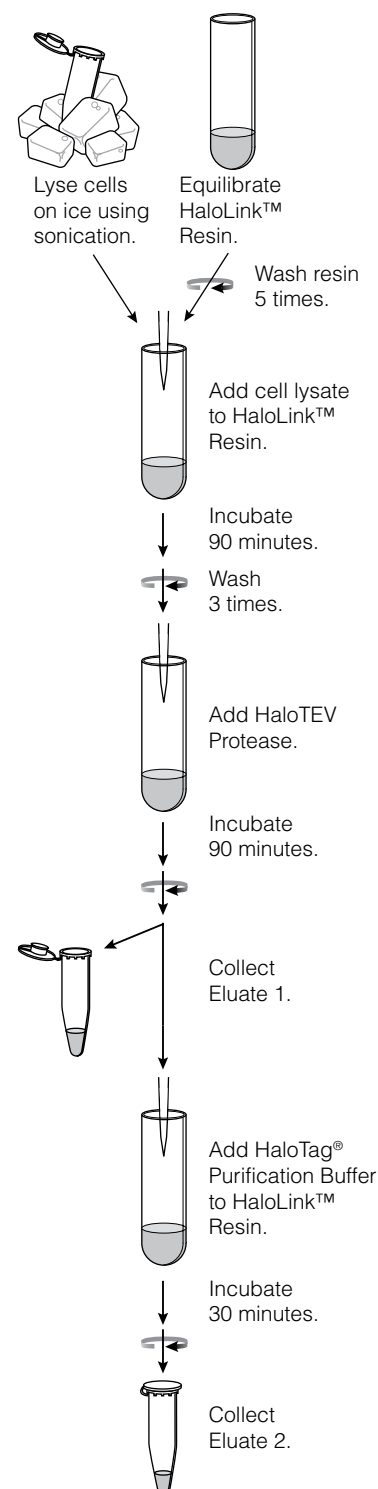
11. Wash the resin three times:
 - a. Add 5ml of HaloTag[®] Purification Buffer; mix at room temperature for 10 minutes.
 - b. Centrifuge at 1,500 \times g for 5 minutes; discard the supernatant.

Cleave

12. Combine 9 μ l of HaloTEV Protease with 291 μ l of HaloTag[®] Purification Buffer.
13. Add the cleavage solution to the resin; incubate at room temperature (22–25°C) for 90 minutes with constant mixing.

Elute

14. Centrifuge at 1,500 \times g for 5 minutes; collect the supernatant (Elution 1).
15. Add 300 μ l HaloTag[®] Purification Buffer to the resin; mix for 30 minutes at room temperature.
16. Transfer the resin into the spin column; centrifuge at 10,000 \times g for 15 seconds; collect Elution 2.
17. Centrifuge Elution 1 and Elution 2 at 10,000 \times g for 1 minute, and transfer to clean tubes.



96061MA

HaloTag[®] Mammalian Protein Detection and Purification Systems

Quick Protocol

Instructions for Use of Products **G6790** and **G6795**

HaloTag[®] Fusion Protein Detection

Fluorescent Labeling

Fluorescent labeling of HaloTag[®] fusion protein with the HaloTag[®] TMRDirect[™] Ligand provides a rapid and convenient method to monitor protein expression and follow the purification efficiency.

1. Dilute the HaloTag[®] TMRDirect[™] Ligand stock solution (100 μ M) twofold in DMSO to make a 50 μ M working solution. Store protected from light, at -20°C .

Note: Alternatively, the stock solution can be prepared in PBS, but cannot be stored.

2. Combine 10 μ l of lysate containing the HaloTag[®] fusion protein with 19 μ l of HaloTag[®] Protein Purification Buffer and 1 μ l of 50 μ M HaloTag[®] TMRDirect[™] Ligand.

Note: The equivalent amount of unbound fraction can be added in place of the lysate.

3. Incubate at room temperature for 15 minutes protected from light.
4. Add 10 μ l of 4X SDS gel loading buffer and heat at 70°C for 3 minutes.
5. Load 10 μ l onto an SDS-polyacrylamide gel.
6. Following electrophoresis, scan the gel on a fluorescence imager using settings appropriate for the HaloTag[®] TMRDirect[™] Ligand (555nm excitation, 585nm emission), and quantitate band intensities.

For further information regarding HaloTag[®] labeling, refer to the *HaloTag[®] Technology: Focus on Fluorescent Imaging with DMSO-Soluble Ligands Technical Manual #TM260* or visit:

www.promega.com/protocols

