
His and GST Purification

***E. coli* Control Lysate**

**Catalog Number
620-0233**

BIO-RAD

Quantity: ~ 30 mg of total protein; ~ 10 mg of target protein.

Description: The His and GST Purification *E. coli* control lysate contains a lyophilized extract prepared from an *E. coli* derivative host. The control lysate contains ~10 mg of a 51 kD dual-tagged target protein. The target protein can be purified using Profinia native IMAC or GST methods. The target protein cannot be purified using denaturing IMAC methods.

Storage: Store at 4°C.

Instructions for Use: Rehydrate the entire vial with 12 ml of 1x native IMAC or GST lysis buffer (catalog #620-0205 or 620-0213) and swirl to mix. Do not shake. Let the vial completely rehydrate for 5 min at room temperature. The rehydrated vial provides enough material for two 1 ml affinity purifications. The prepared lysate can be stored at -20°C for up to 6 months and thawed once prior to use. Do not subject the con-

trol lysate or unused solutions to multiple freeze/thaw cycles.

Shelf Life: The unopened control lysate is stable for at least one year from date of receipt. Refer to the specific expiration date on the product label.

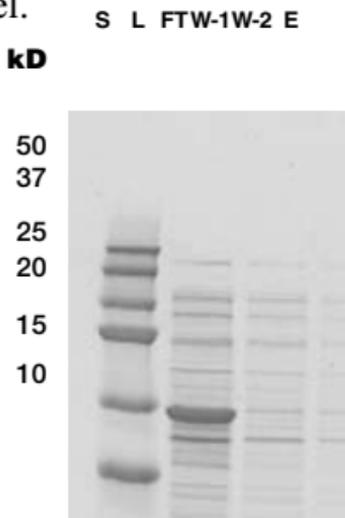


Fig. 1. 51-D IMAC purification. Lanes left to right, Precision Plus Protein™ standards (S); load (L); flowthrough (FT); wash 1 (W-1); wash 2 (W-2); eluted protein (E).

