



# Automated MagneHis™ Protein Purification System

## Automated Protocol #EP011

DESCRIPTION OF THE BECKMAN COULTER BIOMEK® 2000 AND BIOMEK® FX, AND TECAN GENESIS® AND FREEDOM EVO™ METHODS FOR PRODUCT V8550. Please visit the web site to verify that you are using the most current version of this Automated Protocol.

- I. Description ..... 1
- II. Product Components ..... 2
- III. Before You Begin ..... 2
- IV. Automated Processing Requirements for the Biomek® FX ..... 3
  - A. Instrument Requirements for the Biomek® FX ..... 3
  - B. Labware Requirements for the Biomek® FX ..... 3
  - C. Initial Deck Layout for the Biomek® FX ..... 4
  - D. Biomek® FX Specific Pre-Run Recommendations ..... 5
- V. Automated Processing Requirements for the Biomek® 2000 ..... 5
  - A. Instrument Requirements for the Biomek® 2000 ..... 5
  - B. Labware Requirements for the Biomek® 2000 ..... 5
  - C. Initial Deck Layout for the Biomek® 2000 ..... 6
  - D. Reagent Dispense Volumes for the Biomek® 2000 ..... 6
- VI. Automated Processing Requirements for the Tecan Genesis® and Freedom EVO™ Workstations ..... 7
  - A. Instrument Requirements for the Tecan Workstations ..... 7
  - B. Labware Requirements for the Tecan Workstations ..... 7
  - C. Initial Deck Layout for the Tecan Workstations ..... 8
  - D. Tecan Workstation Specific Pre-Run Recommendations ..... 9
- VII. Description of Automated MagneHis™ Protein Purification System ..... 9
- VIII. General Guidelines for Adaptation to Alternative Robotic Platforms ..... 9

### I. Description

This document describes automation of the MagneHis™ Protein Purification System(a,b,c). Specific instructions are provided for the Beckman Coulter Biomek® FX and Biomek® 2000, and Tecan Genesis® and Freedom EVO™ automated liquid-handling workstations. Validated methods for these liquid-handling workstations are available at: [www.promega.com/automethods/](http://www.promega.com/automethods/). General automation guidelines are provided for adaptation to other liquid handling platforms. For troubleshooting chemistry issues please refer to the *MagneHis™ Protein Purification System Technical Manual #TM060*.

**Note:** All Promega Technical Bulletins are available at: [www.promega.com/tbs/](http://www.promega.com/tbs/)

 **Do not freeze**  
the MagneHis™  
Ni-Particles.

## II. Product Components

Product	Size	Cat.#
MagneHis™ Protein Purification System	10ml	V8550

Includes:

- 250ml MagneHis™ Binding/Wash Buffer
- 50ml MagneHis™ Elution Buffer
- 60ml FastBreak™ Cell Lysis Reagent, 10X
- 10ml MagneHis™ Ni-Particles
- 2 vials DNase I (lyophilized)
- 1 Protocol

**Storage Conditions:** Store all MagneHis™ Protein Purification System components at 4°C. **Do not freeze the MagneHis™ Ni-Particles.** The DNase I (lyophilized) may be stored at room temperature. Upon resuspension in water, store the DNase I in aliquots at –20°C. For convenience, the resuspended DNase I may be stored at 4°C for up to one week. **FastBreak™ Cell Lysis Reagent** may form a precipitate at low temperature. If this occurs, warm the reagent to room temperature before use.

## III. Before You Begin

### Materials to Be Supplied by the User

- Tabletop centrifuge capable of 1,500 x g.
- Deep Well MagnaBot® 96 Magnetic Separation Device (Promega Cat.# V3031)
- 1/4 inch Foam Spacer (Cat.# Z3301)
- 2.2ml deep-well, square-well culture plates (Promega Cat.# V6781)
- 1.2ml deep-well, square-well culture plates (Promega Cat.# V6771)
- U-bottom multiwell plates (Promega Cat.# A9161)

### For Biomek® FX:

- Pyramid-bottom reservoir plates (Promega Cat.# V6801)
- Beckman Coulter Orbital Shaker ALP

### For Biomek® 2000:

- Twinivair low-volume, quarter-vertical reservoir (Acme-Automation Cat.# C5001)
- Variomag Teleshake (3mm orbit) (Promega Cat.# V6751)
- Shaker Integration Plate for Biomek 2000 (Promega Cat.# V3691)

### For Tecan Workstations:

- Tecan Te-Shake Orbital shaker (3mm orbit)

**Sample Preparation Before Automated Processing:** If the bacterial cell pellets are frozen, thaw at room temperature for 10–15 minutes before use.

For cell pellets, dilute the supplied FastBreak™ Cell Lysis Reagent to 1X concentration with water before use. For cell cultures, the supplied 10X FastBreak™ Reagent is added directly to the cell cultures at a 1:10 ratio (i.e., 100µl 10X Fastbreak™ Reagent to 900µl culture volume).

**IV. Automated Processing Requirements for the Beckman Biomek® FX**
**A. Instrument Requirements for the Biomek® FX**

<b>Part Description</b>	<b>Quantity</b>	<b>Beckman Coulter Part Number</b>
Minimum: Biomek® FX software version 2.5		Contact Beckman Coulter
96-channel POD	1	Contact Beckman Coulter
Labware positions for 1 POD	11	Contact Beckman Coulter
Tip loader ALP	1	719356
Orbital shaker ALP (optional)	1	Contact Beckman

**B. Labware Requirements for the Biomek® FX**

<b>Part Description</b>	<b>Quantity</b>	<b>Part Number</b>
Deep Well MagnaBot® 96 Magnetic Separation Device	1	Promega Cat.# V3031
2.2ml deep-well plate or comparable	2	Promega Cat.# V6781
1/4 inch Foam Spacer	1	Promega Cat.# Z3301
1.2ml deep-well plate (or comparable)	1	Promega Cat.# V6771
Pyramid-bottom 96-well reservoirs (or comparable)	3	Promega Cat.# V6801
Polystyrene U-bottom multiwell plates (or comparable)	2	Promega Cat.# A9161
200 tips-barrier (rack)	3	Beckman Cat.# 717253

### C. Initial Deck Layout for the Biomek® FX

This is an example of the MagneHis™ Protein Purification System deck layout on a Biomek® FX. Your specific deck layout may be different depending on your Biomek® FX configuration.

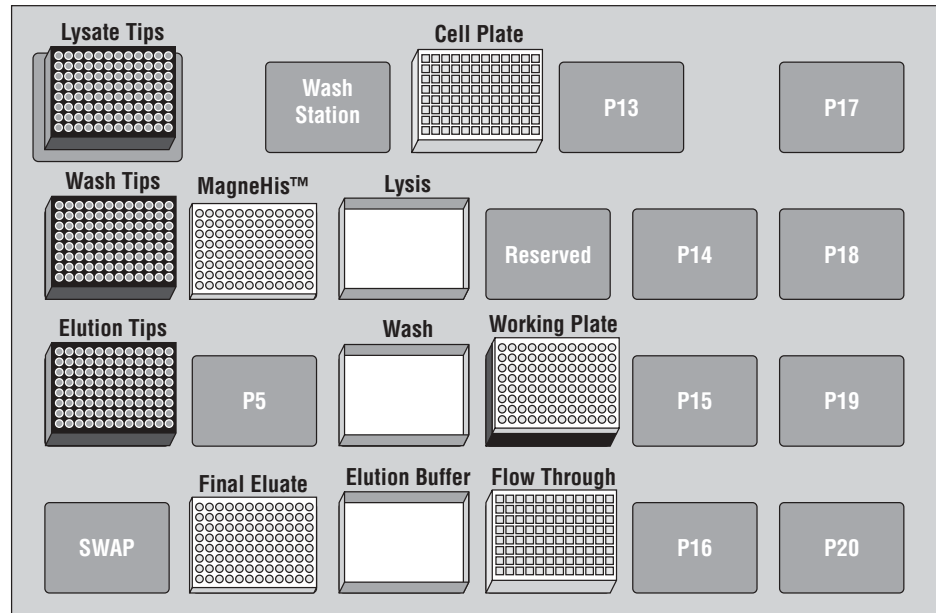


Figure 1. Initial deck configuration of the Beckman Biomek® FX.

ALP Name	Equipment
Tip Loader	200µl ART® Biomek® FX Tips
P1	200µl ART® Biomek® FX Tips
P2	200µl ART® Biomek® FX Tips
P3	Empty: Tip box swap spot
WS1	Tip Wash Station (96 channel; not necessary for method)
P4	96-well, round-bottom plate containing 45µl of MagneHis™ Resin/well (labeled "MagneHis")
P5	Empty
P6	Empty 96-well round-bottom plate (labeled "Final Eluate")
P7	Pyramid-bottom reservoir plate containing 25ml FastBreak™ Cell Lysis Reagent (10X concentration for cultures; 1X concentration for cell pellets) + 25µl DNase I, if desired (labeled "Lysis")
P8	Pyramid-bottom reservoir plate containing 50ml MagneHis™ Binding/Wash Buffer (labeled "Wash")
P9	Pyramid-bottom reservoir plate containing 25ml MagneHis™ Elution Buffer (labeled "Elution")
P10	Empty. Reserved for plate movement
P11	Empty 1.2ml 96-well, round-bottom plate (labeled "Working Plate") on a Deep Well MagnaBot® 96 Magnetic Separation Device with 1/4 inch Spacer
P12	Empty 2.2ml deep-well square-well plate to collect flowthrough (labeled Flow-thru")
Orbital 1	2.2ml deep-well square-well culture plate containing 900µl cell culture in suspension or cell pellets (labeled "Cell Plate")

#### D. Biomek® FX Specific Pre-Run Recommendations

The Biomek FX® automated platform allows users the flexibility to configure the robot's deck according to need. Because of this flexibility in deck configuration, it is likely that the deck used for writing a Biomek® FX method will differ from an end-user's deck. Therefore, it will be generally necessary to map an imported method onto an end-user's deck configuration. To do this, follow the instructions provided in the document **Deck Mapping on the Biomek® FX** ([www.promega.com/automethods/beckman/biomekfx/default.asp](http://www.promega.com/automethods/beckman/biomekfx/default.asp)).

#### V. Automated Processing Requirements for the Beckman Biomek® 2000

##### A. Instrument Requirements for the Biomek® 2000

Part Description	Quantity	Beckman Coulter Part Number
Biomek® 2000 Workstation, 50/60 Hz, 100-120V	1	609000
Biomek® 2000 controller NT	1	609875
BioWorks 3.2 for Beckman Coulter computer	1	609983
MP200 pipetting tool	1	609025
Gripper tool system for Biomek® 2000	1	609001
Tip rack holder	2	609121
Gray labware holder	6	609120
Reservoir holder	1	372795
Quarter vertical reservoir	1	372788
Quarter single reservoir	2	372790

##### B. Labware Requirements for the Biomek® 2000

Part Description	Quantity	Ordering Information
Deep Well MagnaBot® 96		
Magnetic Separation Device	1	Promega Cat.# V3031
1/4 inch Foam Spacer	1	Promega Cat.# Z3301
2.2ml deep-well plate or comparable	2	Promega Cat.# V6781
1.2ml deep-well plate	1	Promega Cat.# V6771
Polystyrene U-bottom multiwell plates or comparable	1	Promega Cat.# A9161
Biomek® 2000 standard 250 tips (rack)	2	Beckman Cat.# 717253
Twinivoir low volume quarter vertical reservoir	1	Acme-Automation Cat.# C5001

### C. Initial Deck Layout for the Biomek® 2000

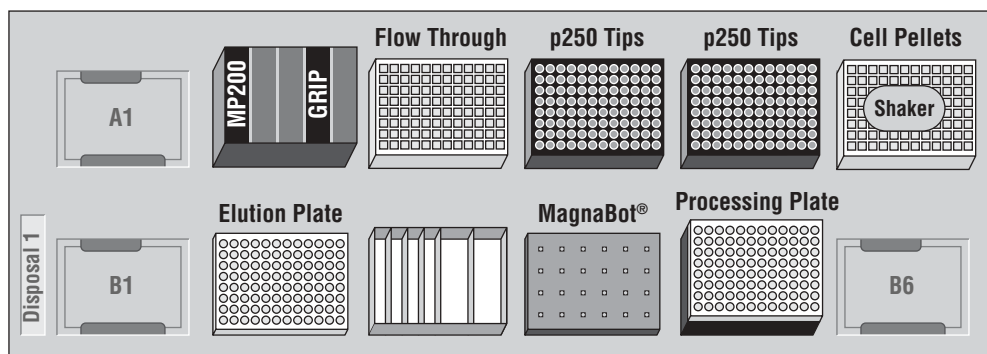


Figure 2. Beckman Biomek® 2000 initial deck configuration.

Position	Equipment
A1	Empty (not used in method)
A2	Tool rack containing MP200 and Gripper tools
A3	Labware holder, 2.2ml deep-well, square-well plate for flowthrough (labeled "Flow Thru")
A4	Tip rack holder, P250 tips
A5	Tip rack holder, P250 tips
A6	Teleshake. 2.2ml deep-well, square-well plate containing samples (labeled "Culture Plate" or "Cell Pellets")
B1	Empty (not used in method)
B2	Labware holder, empty 96-well U-bottom plate (labeled "Elution Plate")
B3	Labware holder, reservoir holder, 1 quarter vertical reservoir, 1 Twinivoir reservoir, and 2 quarter single reservoirs
B4	Labware holder, Deep Well MagnaBot® 96 Magnetic Separation device with 1/4 inch Foam Spacer
B5	Labware holder, empty 1.2ml deep-well U-bottom plate (labeled "Processing Plate")
B6	Labware holder

### D. Reagent Dispense Volumes for the Biomek® 2000

Prior to beginning run, the following MagneHis™ Protein Purification System reagents need to be dispensed appropriately on the deck of the Biomek® 2000:

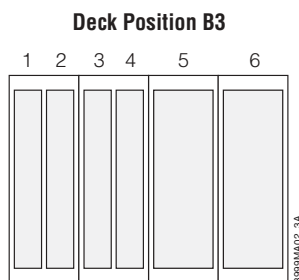


Figure 3. Reagent dispense volumes at Deck Position B3.

- 11ml Elution Buffer
- 11ml 10X FastBreak™ Cell Lysis Reagent (for cell cultures)  
or 22ml 1X FastBreak™ Cell Lysis Reagent (for cell pellets)  
25µl DNase I may be added to the FastBreak™ Reagent if desired.
- 4ml MagneHis™ Ni-Particles
- Empty
- 45ml Binding/Wash Buffer
- Empty (for waste collection)

## VI. Automated Processing Requirements for the Tecan Genesis® and Freedom EVO™ Workstations

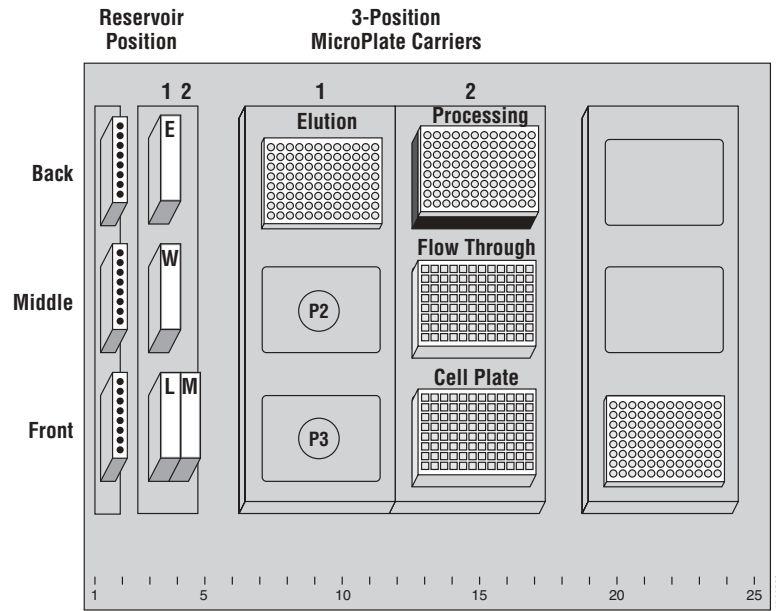
### A. Instrument Requirements for the Tecan Workstations

Part Description	Quantity	Tecan Part Number
Tecan Genesis RSP100 RSP150 or RSP200 or Freedom EVO™ 100, 150 or 200	1	Contact Tecan
Gemini 4.2 software	1	Contact Tecan
1ml syringes	8	Contact Tecan
Standard fixed tips	8	612501
100ml reservoir carriers	2	613020
100ml reservoirs	4	613021
Robotic manipulator arm (RoMa)	1	Contact Tecan
Te-Shake orbital shaker (1 or 2 position)	1	Contact Tecan
3-position microplate carrier	2	Contact Tecan

### B. Labware Requirements for the Tecan Workstations

Part Description	Quantity	Ordering Information
Deep-well MagnaBot® 96		
Magnetic Separation Device	1	Promega Cat.# V3031
1/4 inch Foam Spacer	1	Promega Cat.# Z3301
1.2ml deep-well plate	1	Promega Cat.# V6771
MagnaBot® Adaptor T1	1	Promega Cat.# V8481
2.2ml deep-well plate (or comparable)	2	Promega Cat.# V6781
Polystyrene U-bottom multiwell plate (or comparable)	1	Promega Cat.# A9161

### C. Initial Deck Layout for the Tecan Workstations



**Figure 5. Tecan Genesis® and Freedom EVO™ initial deck configuration.**

Reservoir Position	Column 1	Column 2
Back	15ml MagneHis™ Elution (labeled E)	Empty
Middle	35ml MagneHis™ Binding/Wash (labeled W)	Empty
Front	25ml 1X FastBreak™ Lysis Reagent (for cell pellets) <b>or</b> 15ml 10X FastBreak™ Lysis Reagent (for cell cultures) (labeled L) + 25µl DNase I, if desired	10ml MagneHis™ Resin (labeled M)

Te-Shake Position	Rack on Carrier
Back	Empty
Front	Empty (deck image shows a plate for pipetting to later)

3-Position Microplate Carrier	Rack on Carrier
Position 1	Empty U-bottom multiwell plate (labeled "Elution")
Position 2	Empty
Position 3	Empty

3-Position Microplate Carrier	Rack on Carrier
Position 1	Empty 1.2ml deep-well plate on MagnaBot® T1 Adaptor, Deep Well MagnaBot® 96 Magnetic Separation Device with 1/4 inch Foam Spacer
Position 2	2.2ml deep-well, square-well plate (labeled "Flow-through")
Position 3	2.2ml deep-well plate containing either cell pellets or cell culture.



#### D. Tecan Workstation Specific Pre-Run Recommendations

Due to differences in instrument reference positions, carriers, plate types and Te-Shake configurations, it is recommended to check and adjust the following settings prior to running the method:

- a. X,Y,Z coordinates for each plate and carrier type used in the method
- b. RoMa vectors
- c. Te-Shake: Orbit = 3mm; Counterweight = none

#### VII. Description of Automated MagneHis™ Protein Purification System

This overview describes general liquid handling steps required for the automated MagneHis™ Protein Purification System and can be adapted to a variety of automated liquid-handling robots. For additional information about adaptation to liquid handling robots other than those referenced above, please see Section VIII, General Guidelines for Adaptation to Alternative Robotic Platforms.

1. **MagneHis™ Particle Dispense.** MagneHis™ Ni-Particles are mixed in the reservoir, then 30µl aliquots are dispensed into a 1.2ml deep-well plate.
2. **Cell Pellet Resuspension/Lysis.** Cell pellets contained in the deep-well 96-square-well culture plate are resuspended with FastBreak™ Cell Lysis Reagent by a combination of mixing and shaking for 5 minutes. If processing cell pellets, 200µl 1X FastBreak™ Reagent is used. If processing cell cultures, 100µl 10X FastBreak™ Reagent is used.  
**Note:** to overcome sticky or 'slimy' cell lysates add 1µl/well of resuspended DNase I.
3. **Bind Protein to MagneHis™ Ni-Particles.** The lysate is added to the MagneHis™ Ni-Particles to initiate binding of His-tagged proteins. The lysate is mixed well with the particles by shaking for 1 minute. After shaking, the plate is placed onto the Deep Well MagnaBot® Device to capture the particles. The supernatant, containing unbound materials, can be saved for analysis or discarded to waste.
4. **Washes.** Three washes are performed with the MagneHis™ Binding/Wash Buffer to remove residual impurities from the sample. During each wash, 100–150µl of MagneHis™ Binding/Wash Buffer is added to the MagneHis™ Ni-Particles. The particles are mixed well by shaking for 1 minute. After shaking, the particles are captured on the Deep Well MagnaBot® Device, and the supernatant is discarded to waste. This process is repeated 2 more times.
5. **Elution.** The samples are eluted from the particles by adding 100µl of MagneHis™ Elution Buffer and mixing well for 1 minute on the shaker. After shaking, the particles are captured on the Deep Well MagnaBot® Device, and the supernatant is saved in a collection plate for analysis.

#### VIII. General Guidelines for Adaptation to Alternative Robotic Platforms

The MagneHis™ Ni-Particles settle over time. It is recommended to thoroughly mix the MagneHis™ Ni-Particles on the automated platform prior to dispensing into samples. Resuspension of the MagneHis™ Ni-Particles can be accomplished by thorough mixing or shaking.



(a) Patent Pending.

(b) Certain applications of this product are covered by patents issued and applicable in certain countries. Because purchase of this product does not include a license to perform any patented application, users of this product may be required to obtain a patent license depending upon the particular application and country in which the product is used.

(c) This product is licensed for use under U.S. Pat. No. 6,174,704.

© 2003, 2005 Promega Corporation. All Rights Reserved.

MagnaBot is a trademark of Promega Corporation and is registered with the U.S. Patent and Trademark Office. MagneHis is a trademark of Promega Corporation.

Biomek is a registered trademark of Beckman Instruments, Inc. BioWorks is a trademark of Beckman Coulter, Inc.

Genesis is a registered trademark of Tecan AG Corporation.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

<b>Promega Corporation</b>	
2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Fax	608-277-2516
Internet	<a href="http://www.promega.com">www.promega.com</a>