

# Assurance® GDS Pickpen® II

Part No: 73091 (8 magnet), 73090 (1 magnet)

## General Description

The Pickpen® II is designed to transfer magnetic particles when working with BioControl's Assurance® GDS pathogen detection kits.

## Procedure

### A. Loading Disposable Tips

With magnets withdrawn into the carrier sleeves, position the Pickpen® II over Pickpen® II tips (Part No: 73085 or 73087) and push down firmly to seat the tips over the carrier sleeves.

### B. Collection of Magnetic Particles

- Check that the tips are firmly in place by depressing the red lower button to the first stop (Pickpen® – 8M) or moving the lower lever down (Pickpen® – 1M) to extend the magnets. Depressing the black upper button or pushing the lower lever up will retract the magnets.
- With the magnets locked in the extended position, dip the tips into the suspension containing the magnetic particles. Collect particles by stirring gently for the specified amount of time while continually moving up and down from the surface to the bottom of the well.

### C. Transfer of Magnetic Particles

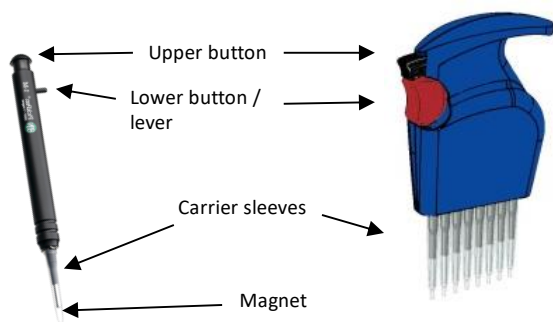
- Once the particles are collected, withdraw the tips from the solution. Tap the Pickpen® II tips against inside of sample wells to remove and residual liquid. The particles will remain attached to the ends of the tips due to the magnetic attraction.
- If necessary, transfer particles to a suitable wash solution to further remove any residual materials.
- Transfer particles to a suitable buffer and depress the black upper button or move the lower lever up to retract the magnets. Gently stir tips in the buffer solution to completely release the particles.

**Note:** To prevent the magnets from not fully retracting. Do not allow fingers to rest on the red button while depressing the black button.

### D. Ejecting Tips

After completing transfer of magnetic particles, the tips can be ejected into a suitable biohazard container by depressing the red button to the second stop (Pickpen® – 8M) or depressing the black upper button (Pickpen® – 1M).

## Pickpen® II – 1M Pickpen® II – 8M



The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

**MERCK**

## Storage

Exposure to temperatures above 80 °C can demagnetize the magnetic tips.

The Pickpen® does not require any calibration and should not be disassembled.

## Cleaning

After use, clean the Pickpen® by lightly spraying a paper towel with 70% isopropyl alcohol and gently wiping the outside of the device. Avoid introducing any liquid into the openings at the end of the carrier sleeves. Exposure of the magnets or the inside of the carrier sleeves to moisture may result in corrosion, preventing the device from operating properly. **Do not spray any solutions directly on the Pickpen®. Do not autoclave.**

If contamination is suspected, moisten paper towel with bleach solution and wipe Pickpen® surface with 10% bleach solution. Allow bleach solution to remain on surfaces for a minimum of 15 min before wiping clean with 70% isopropyl alcohol solution.

To prepare 10% bleach solution add 10 mL of commercially available bleach containing at least 5% sodium hypochlorite to 90 mL of deionized water. The minimum final concentration of sodium hypochlorite in the bleach solution should be 0.5%. The bleach solution is stable for 7 days from preparation. To prepare 70% isopropyl alcohol solution add 70 mL of pure isopropyl alcohol to 30 mL of deionized water or buy commercially available 70% isopropyl alcohol.

## Manufacturing Entity

BioControl Systems, Inc, 12822 SE 32nd St, Bellevue, WA 98005, USA.

BioControl Systems, Inc is an affiliate of Merck KGaA, Darmstadt, Germany.

Merck KGaA  
Frankfurter Strasse 250  
64293 Darmstadt  
Germany  
[Sigmaaldrich.com/GDS](https://Sigmaaldrich.com/GDS)

