

User Guide

Positive Control tool

(Cat. No.: POSICONT1)



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

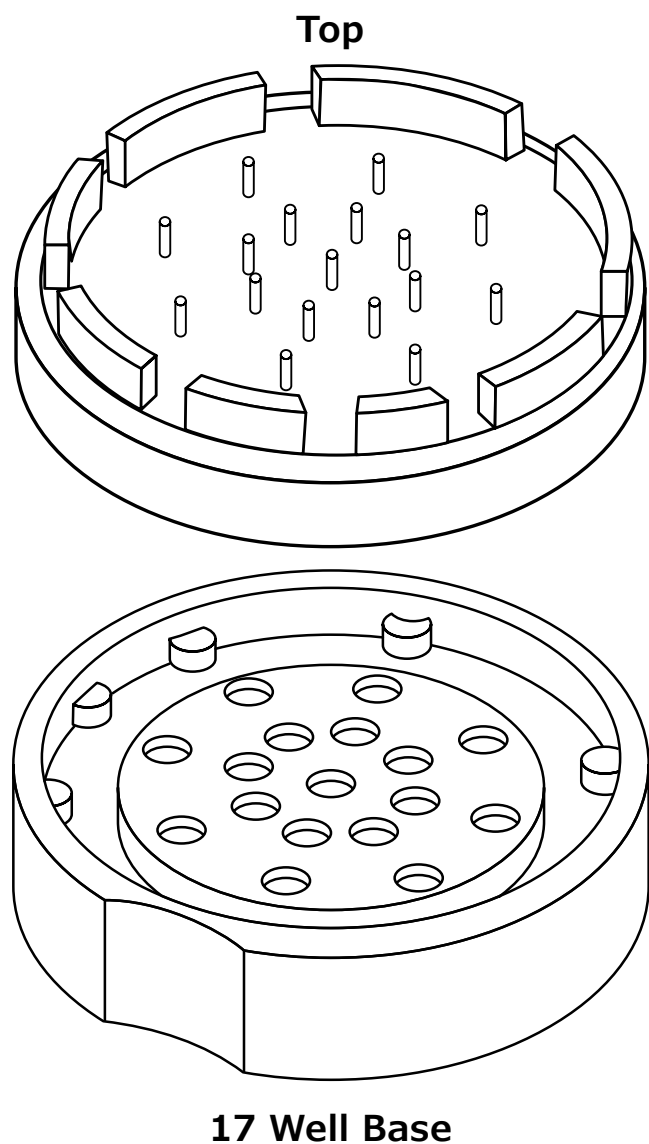
Introduction

The Positive Control Tool is used to verify the functionality of the Milliflex® Rapid System 2.0 AutoSpray Station and the Milliflex® Rapid System 2.0 Detection Tower.

Components

The Positive Control Tool (POSICONT1) consists of:

- Top
- 17 well base



Procedure

Perform the verification of the Milliflex® Rapid AutoSpray Station and the Detection Tower's functionality within a laminar flow hood or a microbial safety cabinet while wearing gloves.

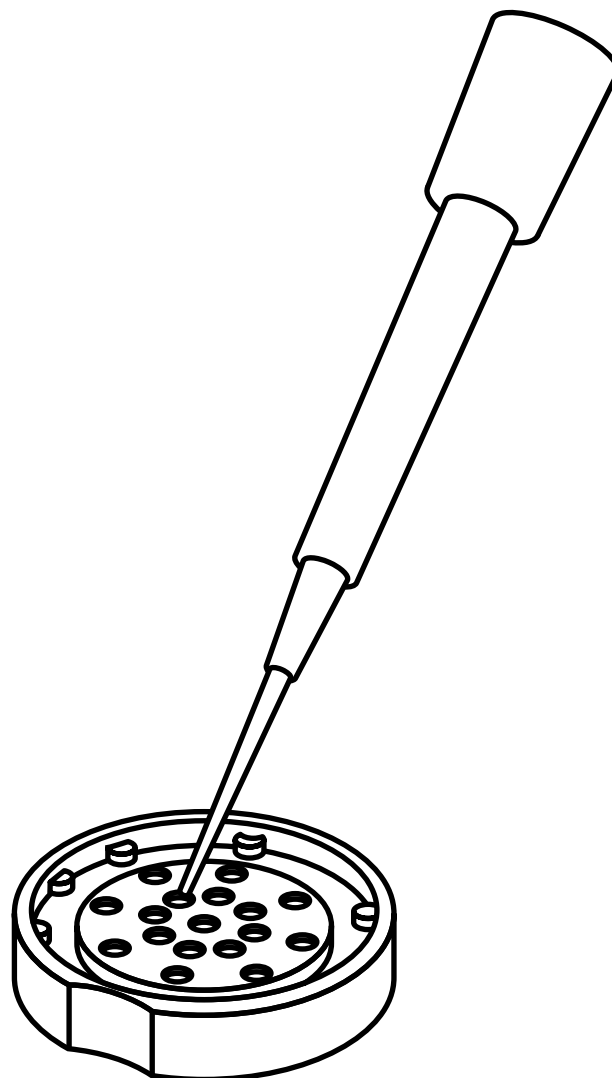
1. Remove the top of the tool.

Note: Ensure that the top does not contact an unclean surface

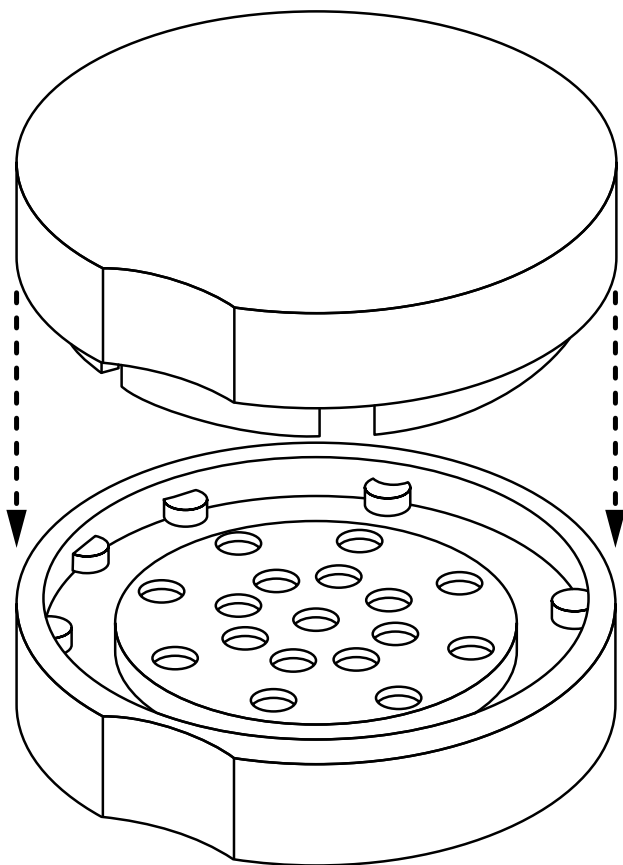
Fill the well(s) of the Positive Control Tool base with an ATP solution as needed. Place 40 µL in each of the 17 wells.

- 2.

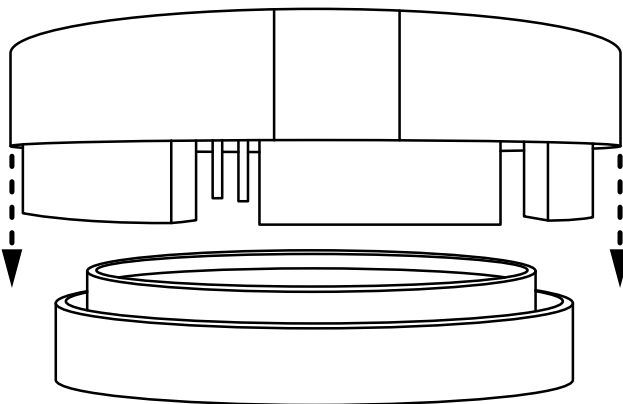
Solution	Concentration
ATP	~10 µg/mL



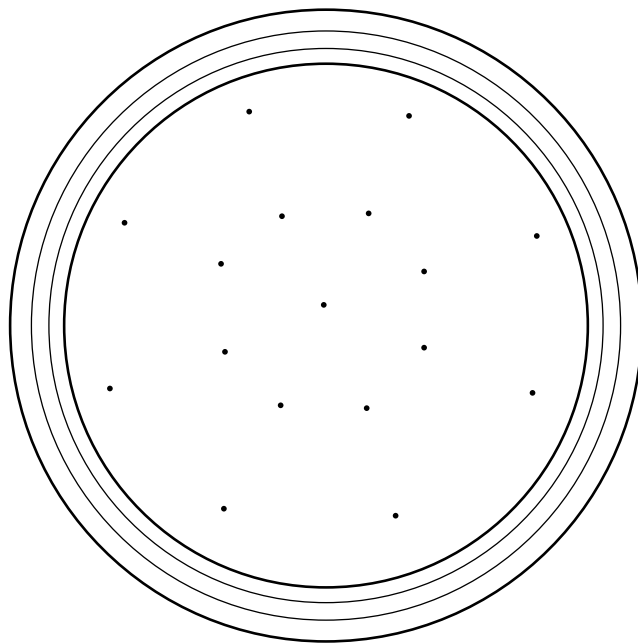
3. To immerse the tips in the solution, align the indent on the top with the indent in the base.



4. Remove the top from the base. Gently install the top of the tool onto the Milliflex® membrane to spot the membrane surface. Ensure that the top maintains contact with the membrane for two seconds.



5. Lift the top straight up to remove it from the membrane. Verify that one or more spots are visible on the membrane.



6. Once the membrane has dried for one minute, spray reagent on it with the AutoSpray Station (refer to the Milliflex® Rapid System 2.0 AutoSpray Station and Milliflex® Rapid System 2.0 Detection Tower User Guides). Subsequently, analyze the membrane using the Milliflex® Rapid System (see the Milliflex® Rapid System 2.0 AutoSpray Station user guide UG11631 and Detection Tower user guide UG11577). The count result must be ≥ 1 count.

Cleaning and Decontamination

The Positive Control Tool is made from aluminum and must be cleaned after each use of ATP solution.

1. Empty the well(s).
2. Rinse with purified water.
3. Autoclave at 121 °C for 30 minutes or at 134 °C for 10 minutes.

Technical Assistance

For more information, please visit SigmaAldrich.com/offices for up to date worldwide contact information.

Manufacturer

Millipore SAS
39 Route Industrielle de la Hardt
67120 Molsheim, France

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

SigmaAldrich.com/milliflex-rapid-2

