

Syringe Filters with Luer Lock Connection Secure Business Sample Processing for Cancer Diagnostic

Customer Challenge

The medical diagnostic manufacturing company needed syringe filtration devices as part of consumables kit for an instrument. They had the following criteria:

- Simplify bench preparation of radio-labelled compounds used in nuclear medicine for cancer diagnostics
- Secure and personalized sample preparation, which includes a series of liquid transfers, fillings and critical manual handlings
- Safer protocol for both patient and clinician

Solution Provided

During the innovation stage, the diagnostic company used Minisart® syringe filters to vent gases in and out of reagentflasks, allowing clean and sterile air to pass through. At the product development stage, they validated a Luer-lock connection as an effective and secure connection to their instrument and checked the syringe filter for air flow rate and pressure.

Sterile and hydrophobic Minisart® PTFE 0.2 µm, diameter 15 mm, was confirmed as the best option to meet their requirements. Inlet was connected to the tubing via a female Luer-lock and the outlet was connected to the instrument with a male Luer-slip connection. Single-blister packing of products ensured sterility.

Once Minisart® was confirmed to fully comply with the manufacturer requirements, the company contacted Sartorius experts to discuss quantities, market expectations, regulations, supply chain, and quality certifications. The diagnostic manufacturer categorized Minisart® syringe filters as non-critical components. The project also involved Legal, Regulatory Affairs and Procurement from both parties.

Regular follow-ups are carried out to optimize supply chain

At a Glance

- Easy-to-plug syringe filters with Luer lock connection
- Gas venting allowing clean and sterile air to pass through
- Patient safety is not compromised
- Download: [Sartorius Minisart® Selection Guide](#)



Case Profile

Company Type::

Instruments and reagents manufacturer - Diagnostic radiotracers

Industry:

Diagnostic

Company Size:

250 Employees

Outcome

The diagnostic manufacturer successfully brought their innovative instrument to the market. Venting filters were included in a ready-to-use set of consumables and reagents that allowed for a compact and flexible configuration of their instrument.