

Minisart® to Secure Sample Processing for Cancer Diagnostic

The Company is dedicated to advancing radiopharmacy solutions for diagnosing and treating significant health issues, including cancer and cardiovascular diseases. They champion nuclear medicine's potential in healthcare, offering user-friendly in-house radiolabeling systems that enable medical staff to prepare, automate, and administer radiopharmaceuticals safely while minimizing radiation exposure.

Testing Challenge

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

The Solution

- Sterile and hydrophobic venting Minisart® PTFE 0.2 µm, diameter 15 mm
- 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 is kitted by the Company for sterility assurance

The Value

- Venting filters included in a ready-to-use set of consumables and reagents allowing clean and sterile air to pass through
- Patient and clinician safety not compromised
- Easy-to-plug with Luer lock connection



For more information,
visit www.sartorius.com

SARTORIUS