

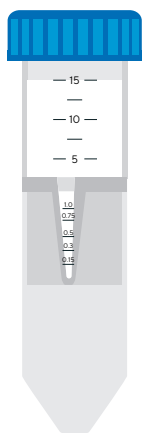
One-Step Milk Fractionation

Efficient Sample Preparation for Residue Analysis

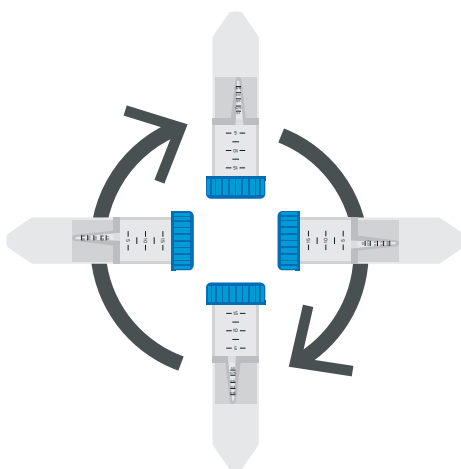
Skip time-consuming methods such as QuEChERS or QuPPe.

Choose ultrafiltration to speed up milk sample preparation for antibiotic, chlorate and perchlorate testing.

Quick and Easy to Use



1 Add raw milk sample



2 Remove fats | proteins



3 LC/MS-ready filtrate



Remove your bottleneck:
Prepare more samples, faster.



Prevent contamination:
Avoid additives to ensure reliable results.



Improve sustainability:
Use fewer reagents and consumables.



Reduce costs:
Make your testing budget go further.

For further information and
to request samples, contact:
office@puremilk.eu

*pure***Milk**
ANALYTICAL

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600



For further information, visit

[sartorius.com](https://www.sartorius.com)