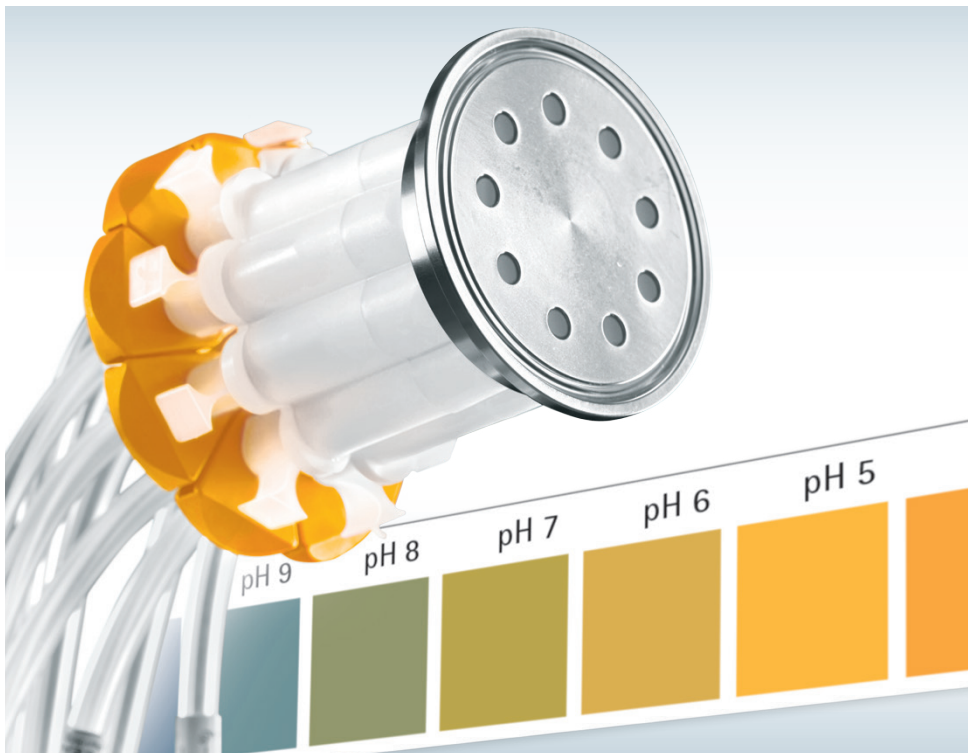




Aseptic Sampling Best Practices: Sample Collection for pH and Gas Analysis



Application
Note

#01

#02

#03

#04

#05

Introduction

Accurate monitoring of dissolved oxygen, carbon dioxide and pH is essential to a successful drug production process.

The most precise results begin with how the sample is collected.

Syringe: Ideal for pH and Gas Analysis

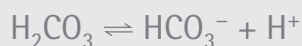
Syringes are an ideal sample collection vessel for pH, DCO_2 and DO_2 analysis. Sartorius consulted with industry leading bioanalyzer manufacturer Nova Biomedical® to detail why.

The Chemistry of pH

1. When carbon dioxide dissolves into water it exists in equilibrium to form carbonic acid:



2. Carbonic acid disassociates to generate a free hydrogen ion which can change a solution's pH:



Time from Sample to Analysis

The pH of a sample will shift as cells within consume O_2 and produce CO_2 . The longer the time between sample collection and analysis the less accurate the analysis test will be.

Samples collected into a syringe can be taken directly to the analyzer for testing, giving the most precise result.

Gas Exchange

De-gassing or off-gassing occurs where the fluid is in contact with a surface or the environment.

Cylinders, like a syringe, maintain a more stable surface area-to-volume ratio. Gas exchange is limited and stabilized.

Materials of Construction

Syringes are constructed of plastic. All plastics will allow some degree of gas transmission into or out of solution. The slower the rate of gas transmission, the more accurate the gas or pH analysis will be.

Syringes are constructed of thick-walled and rigid plastic. Compared to thin-walled flexible plastics gas transmission rates are slowed in a syringe.

Environment

Saturation of O_2 and CO_2 in cell culture can be different from saturation levels in air. Gases in air bind to the plastic surface of the sample collection container prior to use. The bound gases transfer from high to low concentration artificially changing dissolved gas in the sample.

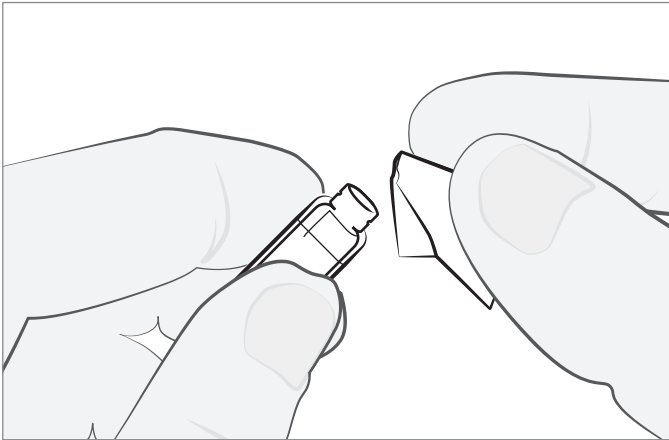
Syringes are not immune to the effect but the more stable surface area-to-volume ratio reduces its impact.

Stable Surface Area to Volume Ratio

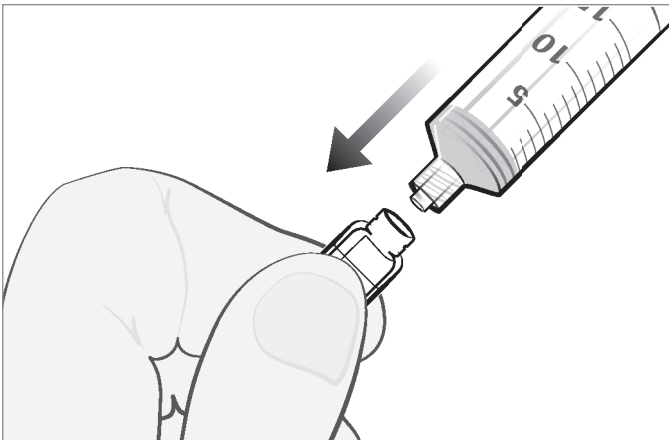
There is more opportunity for gas exchange where the solution is in contact with a surface. Consequently, there is less variability of gas exchange with stable surface area-to-volume ratios.

Sample Volume Collected	10 mL Syringe		50 mL Bag	
	Surface Area (cm ²)	Surface Area to Volume	Surface Area (cm ²)	Surface Area to Volume
2mL	8.82	4.41	424	212.00
5 mL	17.10	3.42	424	84.80
10 mL	30.89	3.09	424	42.40

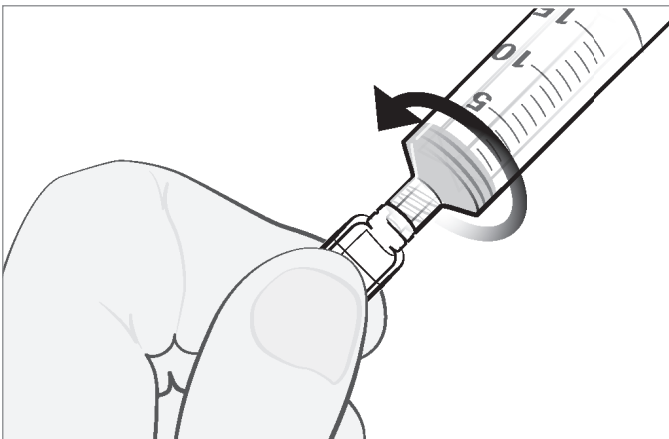
Step-by-Step Instructions for Use



1. Sanitize valve surface



2. Insert fitting straight into valve



3. Turn clockwise to engage

BENCHMARK™ sampling line

- Precise pH, DO₂ & DCO₂ sample analysis
- High frequency sampling from a single sampling line
- Economical solution using conventional equipment and components

Description

BENCHMARK™ features a luer-activated access site which accepts any standard male luer fitting – common to most syringes.

The stem of the male luer opens the silicone valve as it is connected to the access site allowing for unimpeded fluid flow. The valve closes when the male luer is disengaged.

Secure Expansion of Sampling Capacity

The access valve has been validated to maintain a microbial barrier after 140 actuations making it suitable for high-frequency sampling requirements.

Applications

- BENCHMARK™ is ideal for most process-monitoring sampling needs
- Configure BENCHMARK™ on the TAKEONE® aseptic sampling system
- Connect BENCHMARK™ to benchtop bioreactors



Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5 – 7
34302 Guxhagen

Phone +49.5665.407.0
Fax +49.5665.407.2200

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Modcenterstrasse 22
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.647.38.40
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius Stedim RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone +34.913.586.098
Fax +34.913.589.623

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605

Phone +52.55.5562.1102
Fax +52.55.5562.2942

leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Av. Emilio Cavenecia 264 San Isidro
15073 Lima, Perú

Phone +51.1.441 0158
Fax +51.1.422 6100

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Biotech (Beijing) Co. Ltd.
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

Sartorius Stedim Biotech (Beijing) Co. Ltd.
Guangzhou Representative Office
Unit K, Building 23
Huihua Commerce & Trade Building
No. 80 Xianlie Middle Road
Guangzhou 510070

Phone +86.20.37618687 | 37618651
Fax +86.20.37619051

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300
Fax +81.3.4331.4301

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799



▶ www.sartorius-stedim.com