

PHARMACEUTICAL ENGINEERING®



The Official Magazine of ISPE
September–October 2020 | Volume 40, Number 5

The Transition to Digitalization

Validation of Clinical Trial-Related
Systems in Smaller Enterprises

Quality and Regulatory Solutions for
PAT in Continuous Manufacturing

USING CFD MULTIPHASE MODELING
to Predict Bioreactor Performance

SELF-CALIBRATING THERMOMETERS
for Use in Medical Autoclaves



ispe.org/pharmaceutical-engineering

Sartorius Korea Biotech은 세포치료제 개발, 임상연구 및 상업 생산을 위한 **Total Solution** 을 제공합니다.

High-quality serum-free stem cell culture media to support the reproducible generations of quality-assured and clinically acceptable stem cells.



hPSC, MSC, T-CELL, PBMC

Cryopreservation, Recombinant Trypsine, Human Platelet Lysate, Human Serum Albumin, Vitronectin, Attached solution.

- Serum-Free, Xeno-Free, Defined
- Manufactured under cGMP
- DMF (File Master Drug)

Related Products

Regenerative Medicine Bioprocessing



Biostat® RM TX

- Easily scalable expansion Bioreactor with single-use for Cell Therapy



kSep® 400 & 6000S

- Gentle harvesting systems
- Simple, scalable centrifugation



IntelliCyt® iQue3

- Faster and Smarter Flow Cytometry
- Immune cell profiling and QC



New Microsart® ATMP kits

- 3-hour rapid sterility testing transferrable to GMP



2020년 5월, 싸토리우스는 생명 과학 사업 분야에 대한 포트폴리오 확장을 성공적으로 마무리 하였습니다!

인수를 통한 포트폴리오 확장을 통해 바이오 의약품 및 백신과 같은 의약품의 안전하고 효율적인 개발 및 생산에 기여하고, 국내 바이오 산업 전반을 향상 시키는데 이바지 하도록 최선을 다하겠습니다.



Chromatography Systems & Resins

- High selectivity, versatility and efficiency in separating molecules in a mixture



Microcarrier & Cell Culture Media

- Customization, Optimization & Creation



Single-use TFF Systems

- Process Solutions for Ultrafiltration and Diafiltration



Octet Systems

- Superior performance, unrivaled flexibility