



# KOREA BIOPROCESSING EXCELLENCE AWARDS 2020

---

21 - 22 October 2020

**WINNERS HANDBOOK**

BROUGHT TO YOU BY





# KOREA BIOPROCESSING EXCELLENCE AWARDS

The Korea Bioprocessing Excellence Awards seeks to give recognition to exceptional Bioprocessing and Cell & Gene therapy experts, organizations and technologies. Featuring top Bioprocessing and Cell & Gene therapy leaders in the industry, along with the latest advanced technologies and best practices. The Korea Bioprocessing Excellence Awards applauds extraordinary leaders & trend setters of today and inspires innovators of tomorrow.

The Awards Ceremony was held in association with the 9th Biologics Manufacturing Korea and Cell Therapy World Asia 2020 Virtual Conference, which was streamed live on 21 - 22 October 2020.



# CATEGORIES WINNERS

---



Best Biological  
Regenerative  
Medicine Pipeline of  
the year



Bioprocessing  
Excellence in Viral  
Vaccines  
Manufacturing



Bioprocessing  
Excellence in  
Continuous  
Processing  
Implementation



MEDIPOST

Bioprocessing  
Excellence in Cell  
Therapy  
Manufacturing

**SAMSUNG**  
BIOEPIS

Bioprocessing  
Excellence in  
Biosimilars



Bioprocessing  
Excellence in Antiviral  
Therapeutics



Cell therapy  
Bioprocessing  
Excellence of the year

**SARTORIUS**

Best Scalable Cell  
Engineering Platforms  
of the year

**FUJIFILM**  
Value from Innovation

Bioprocessing  
Excellence in Cell  
Culture Optimization  
Technologies in Korea





**MERCK**

Bioprocessing  
Excellence in  
Upstream Processing  
in Korea



**PALL** Biotech

Bioprocessing  
Excellence in  
Downstream  
Processing in Korea



**West**  
By your side  
for a healthier world™

Bioprocessing  
Excellence in Fill-  
Finish Operations



**ThermoFisher**  
SCIENTIFIC  
The world leader in serving science

Bioprocessing  
Excellence in Viral  
Vector Production in  
Korea



charles river

Bioprocessing  
Excellence in Viral  
Clearance and Safety



**MDimune**

Best Cell Therapy  
Technology of the  
Year



**ORGANOID**  
SCIENCES

Best Cell Therapy  
Pipeline of the Year



**Medytox**

Best Cold Chain  
Management in  
Pharma



**SAMSUNG**  
BIOLOGICS

Best Biologics CMO  
of the Year



**avantor**™

Bioprocessing  
Excellence in Single-  
Use Manufacturing in  
Korea



**Envirotainer**®  
The Active Cold Chain

Bioprocessing  
Excellence in Active  
Packaging Solutions  
in Korea



**cytiva**

Best supplier in  
Integrated Downstream  
Bioprocessing in Korea



Innovative  
Bioprocessing  
Technology



Bioprocessing  
Excellence in  
Bioreactors



Bioprocessing  
Excellence in Passive  
Packaging Solutions  
in Korea



Bioprocessing  
Excellence in CDMO  
Services



Bioprocessing  
Excellence in  
Innovative Single-Use  
Solutions in Korea



Bioprocessing  
Excellence in  
Analytical Solutions in  
Korea



Best CMO in Aseptic  
Fill and Finish &  
Packaging in Korea

**Jeonghoon Han**  
Vice President  
Eutilex

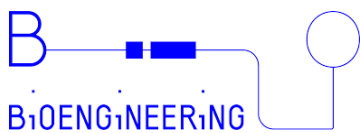


Cell Therapy Lifetime  
Achievement award

**David Bang**  
Managing Partner  
SIG Business Consulting



Bio Logistics Lifetime  
Achievement Award



Bioprocessing  
Excellence in Facility  
Design and  
Engineering Services