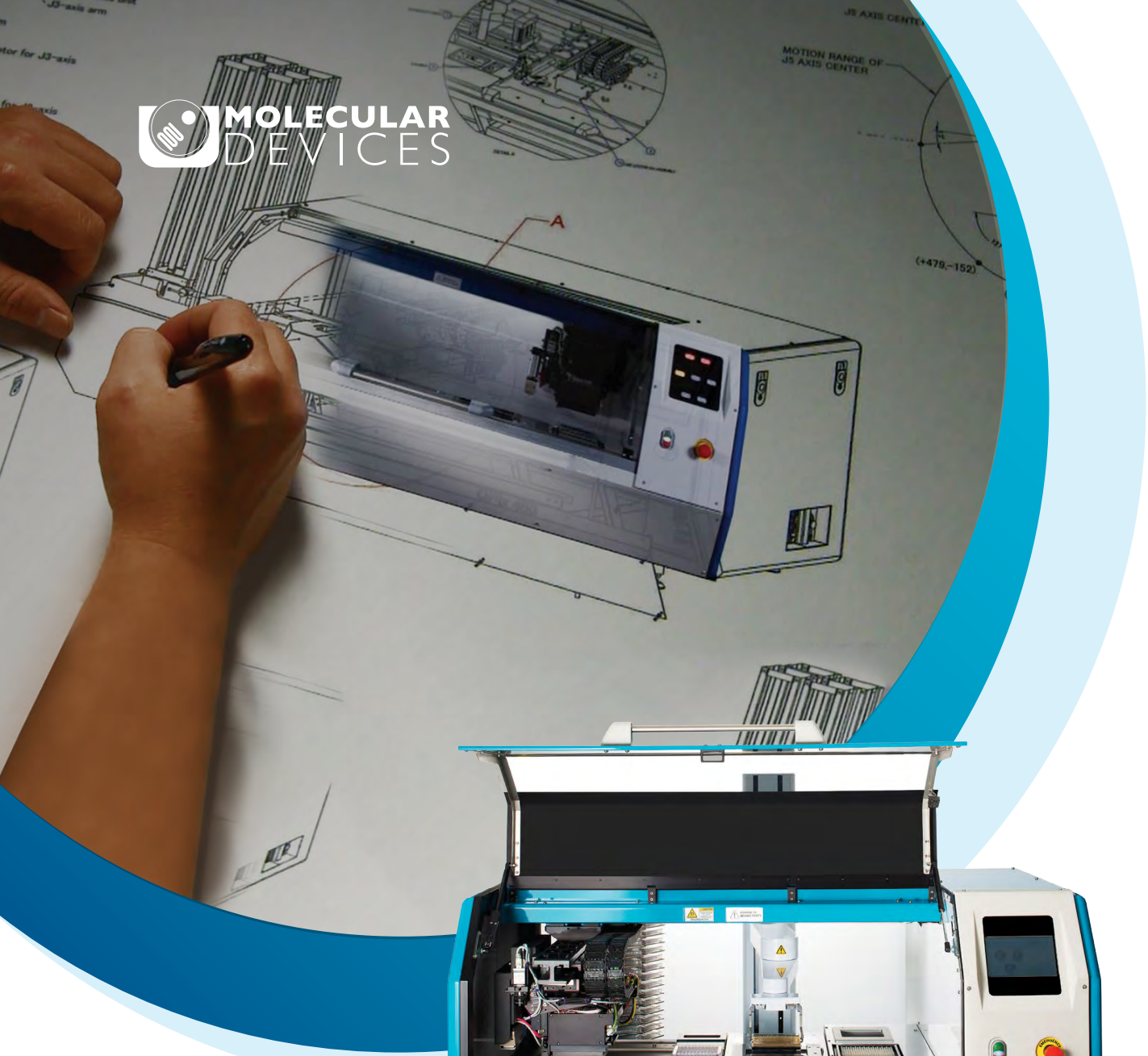


 **MOLECULAR
DEVICES**



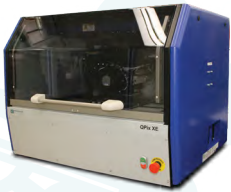
QPix

Microbial Colony Picker

미생물 콜로니 선별 장비

Qpix colony picker

QPix Series 는 Microbiology 관련 연구와 실험 수행을 자동화합니다. 미생물 콜로니를 스크리닝하고 선별하여 Cloning 과정에서의 양성 콜로니 (beta-gal blue colony) 를 자동으로 선별하여 접종하며, 그 외 옵션에 따라 cell plating, spreading, streaking 및 Imaging 기반의 phenotypic selection, cherry picking, library replication & management 등의 기능을 수행하기 때문에, 미생물 관련 워크플로가 진행되는 필요한 대부분의 노동집약적 업무에 효율성과 생산성을 더합니다.



QPix XE system

Compact automated colony picker

컴팩트한 사이즈로 미생물 콜로니 관리의 효율성을 최대화 시킨 모델입니다. 4-12개의 96well plate를 적용할 수 있으며 시간당 최대 1,500 개의 콜로니를 선별합니다.



QPix 420 system

Automated colony picker

Colony screening 배치 규모를 더욱 확장시킨 모델입니다. 12-20개의 96well plate를 적용할 수 있으며, 시간당 최대 3,000개의 콜로니를 선별합니다.



QPix 450/460 systems

Plating to picking

3개의 staker 를 추가하여 최대 210개의 destination plate 를 수용할 수 있습니다. Liquid handling, plating (spreading), streaking 옵션을 추가할 수 있습니다.



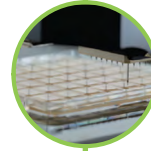
QPix HT system

Flexible, modular, fully automated colony picking and library management system

현재 진행중인 워크플로와 실험 세팅에 맞추어 전반적인 자동화 시스템 구현이 가능합니다. Sample plating, streaking 및 자동화 robotics 를 연결하여 throughput 을 최대화 할 수 있습니다.



Wash Bath + Halogen Lamp
2중 멸균 시스템



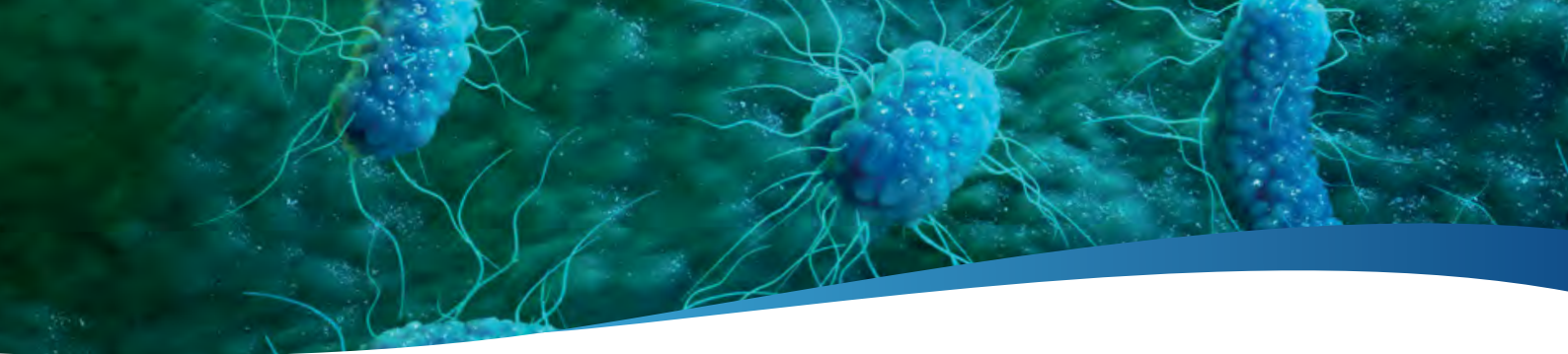
Agar Plate 높이
측정하여 인식



Barcode 리더기
콜로니 추적 가능



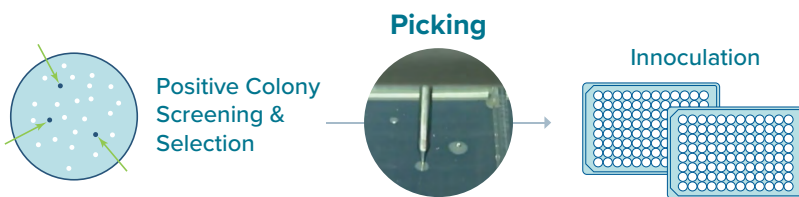
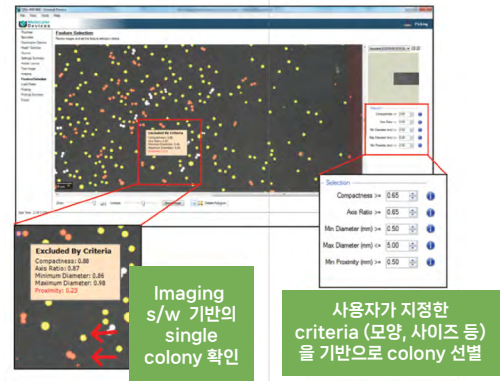
Strain Size에 맞는
Picking Pin 적용 가능



Colony picking

Positive colony selection and inoculation

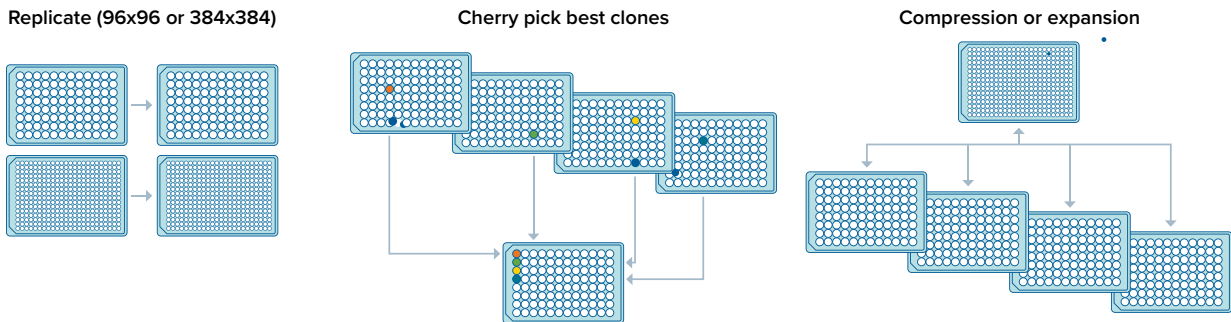
이미징을 기반으로 미생물(박테리아, 진균, 효모, Viral plaque) 군집을 정확하게 스크리닝하고 선별합니다. 한시간에 1500 - 3000 개의 콜로니를 picking 하며 2중 (Halogen + EtOH) 멸균 관리 및 Barcode 를 통한 데이터 관리가 가능합니다.



Library management

Plate replacement and rearrangement

Bacterial Library, Sample Bank 관리에 유용한 Plate Replication, Compression, Expansion 및 Best Clone 의 Cherry Picking 과정을 정확하고 신속하게 자동으로 수행합니다.

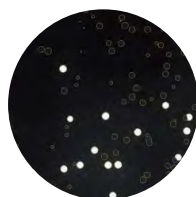


Automated microbial workflow

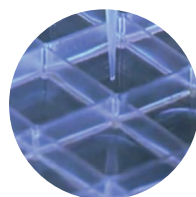
Phenotypic selection, automated cell plating and streaking, plaque and halo detection, etc.



Beta-gal selection
Phenotyping selection
Molecular cloning



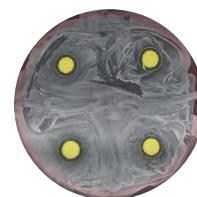
Beta-gal selection
Gene editing
Protein engineering



Protein expression
Seeding/spreading
/streaking



Plaque selection
Phage display



Halo detection
Antibiotics susceptibility

QPix applications

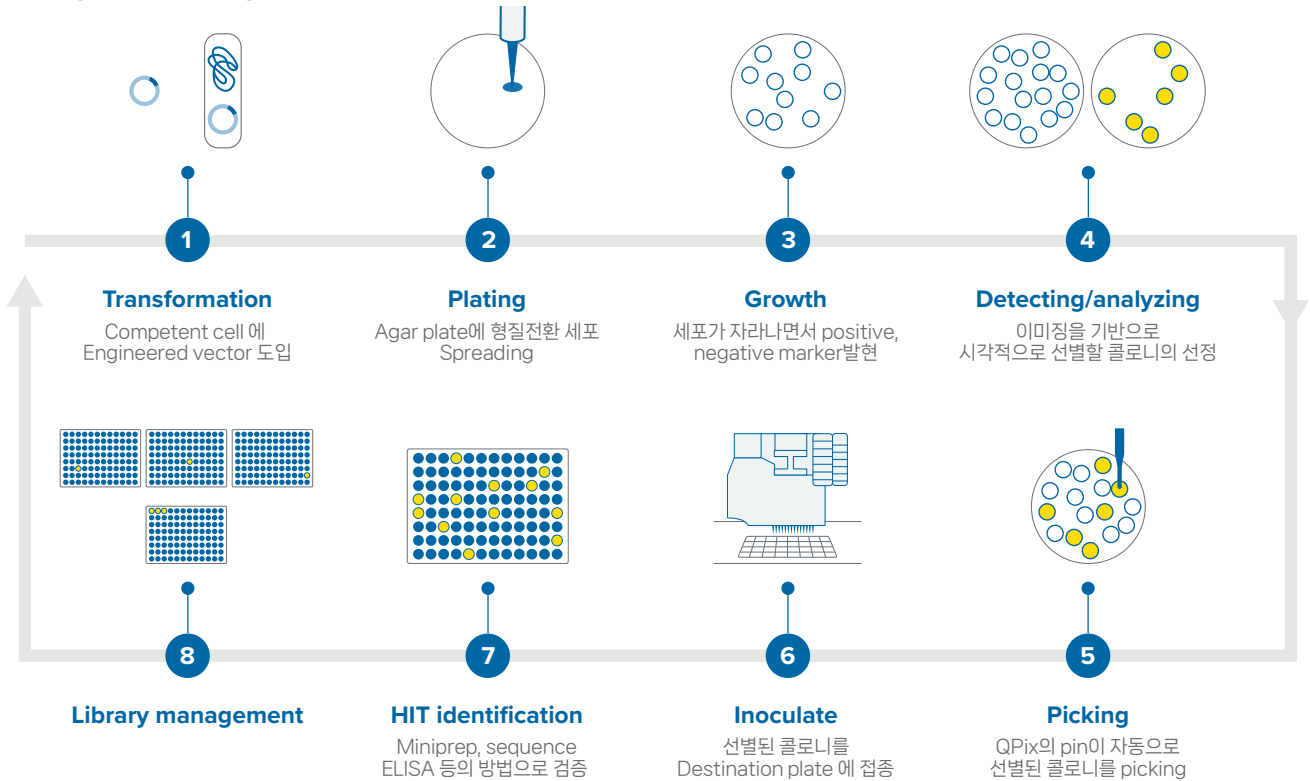
QPix 미생물 콜로니 선별 장비는 미생물을 활용한 실험이나 유전자 재조합이 필요한 다양한 영역에서 사용되고 있습니다.

Protein engineering
Gene editing
Gene cloning
Synthetic biology

Biofuel
Agriculture
Microbiome
Cell line development

Microbial strain engineering
Environmental science
Food and beverage
Phage display

QPix - gene cloning workflow



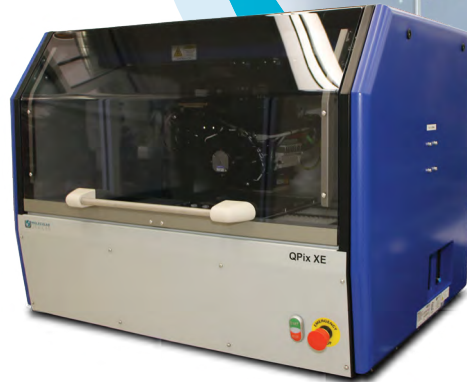
Customer case: Inscripta

1,000-10,000 개의 콜로니 선별작업이 요구되는 digital genome engineering (Onyx) 기술과 QPix 를 효과적으로 접목하여 합성생물학의 워크플로를 혁신적으로 개선한 next generation synthtic biology 솔루션을 제시할 수 있었습니다.



New

QPix XE Automated colony picker



QPix XE는 대학 및 소규모 랩에서 활용 가능한 컴팩트한 사이즈의 자동화 미생물 콜로니 선별 장비입니다. 고급 Imaging 과 Robotic 기능을 탑재하여 박테리아, 효모 및 곰팡이를 포함한 다양한 샘플 유형에서 개별 colony를 정확하고 효율적으로 선별할 수 있습니다. 작은 실험실에서도 공간을 많이 차지하지 않아 효율적이고 이상적인 솔루션입니다.



이미징, Phenotype 기반의
미생물 콜로니 선별 자동화



시간당 최대 1000 - 1500개의
콜로니 자동 처리



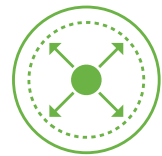
간편한 소프트웨어와 UI 로
쉽고 직관적인 사용법



공간 활용을 최대화 시킨
미니멀한 사이즈



자동 샘플 추적 및
데이터 저장



원하는 Module 을 추가하여
확장 가능성



Colony picking 시
Agar plate 높이 측정 센서 활용



다양한 Labware 및
미생물 종에 적용 가능

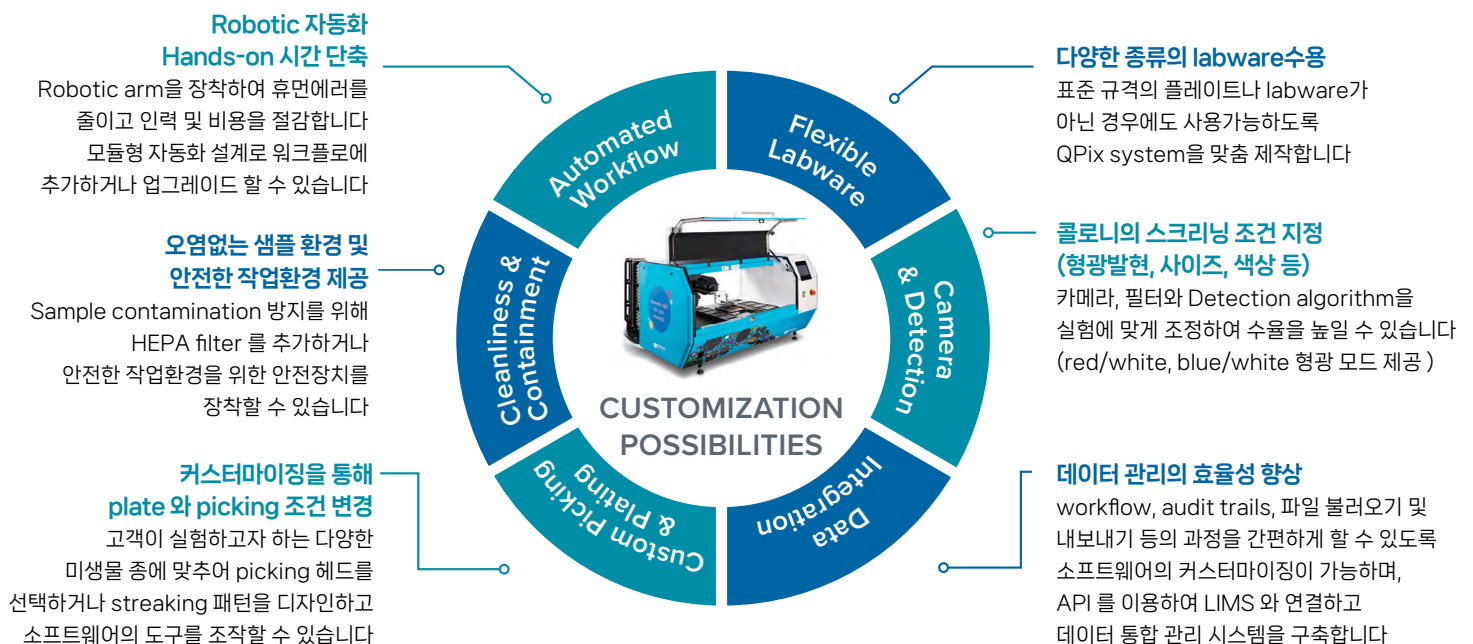


백색광 + 형광(Optional)
이미징 제공



QPix custom microbial colony picker

QPix® Microbial Colony Picker 고객의 실험디자인과 요구사항에 맞추어 커스터마이징 하여 통합 자동화 워크스테이션을 제공합니다.



Custom project 사례

Edinburg Genome Foundry (EGF)

DNA manufacturing 을 위한 자동화 시스템 구축

고객 요청사항:

large DNA design 및 assembly 를 위한 완전 자동화 워크플로를 구현하기 위해 QPix colony picker 와 함께 3개의 robotic arm, 2 liquid handler, incubator, PCR machine 이 들어간 통합 워크스테이션 구축

AWS:

- Robotic arm 이 접근 가능하도록 QPix system 구조 변경
- 8개의 샘플을 동시에 plating 하기 위한 custom picking head 개발
- 원활한 장비 Integration을 위해 open API 및 s/w 지원

결과

맞춤제작된 QPix system은 시간당 3천개의 colony 를 처리하여 효과적인 sample data 관리를 위한 통합된 tracking을 제공하였으며 노동집약적인 업무 단축 및 운영비가 절감됨



Robotic arm의 접근이 가능하도록 Qpix후면 개조



개조된 QPix system (파란 화살표) 부터 plate hotel 까지 접근 가능한 robotic arm

Customer X

Protein engineering 을 위한 라이브러리 스크리닝

고객 요청사항:

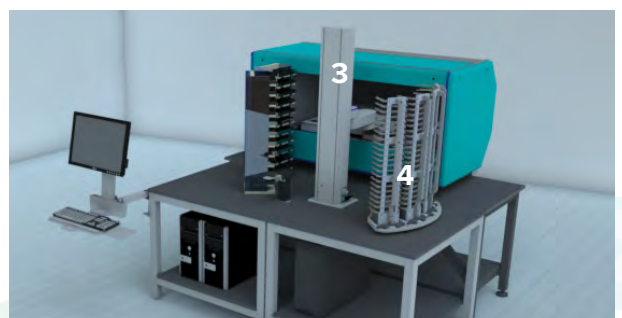
기능성 단백질 생산을 위해 박테리아 라이브러리를 통해 항체와 단백질 스크리닝작업을 필요로 했으며, Colony selection 및 inoculation 과정 및 하루에 30개의 96well plate 및 deep well plate 에 직접적으로 plating 을 처리하는 시스템 필요

AWS:

- Robotic arm이 접근 가능하도록 QPix system 구조 변경
- Green Button Go 소프트웨어, Preciseflex PF3400 robot arm, Ambient plate, QTray storage (88 plate) 와 HEPA 필터가 적용된 에어클리너를 포함하는 통합 워크스테이션 구축

결과

맞춤 제작된 Qpix 시스템으로 시간당 3천 개의 콜로니를 처리하면서 80개 이상의 플레이트를 한번에 처리하여 노동집약적인 업무 단축 및 운영비 절감



(1) Robotic arm의 접근이 가능하도록 개조된 QPix (2) HEPA enclosure (3) PF3400 (4) Plate storage

QPix specification 비교



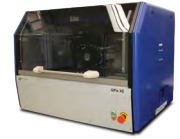
QPix 420



QPix 450/460



QPix HT



QPix XE

	Basic design for automating colony picking with small footprint. Ideal system to replace manual with automated picking and allows for a flexible bed setup and labware use.	From plating to picking – increase throughput with up to 210 destination plates in three stacker lanes. Optional fluidics for plating and streaking allows you to plate and pick samples.	Flexible, modular, and fully automated colony picking and library management system is ready for robotics integration for maximum throughput and walkaway time.	Ideal for small lab spaces with no compromise on efficiency, with advanced imaging and robotic capabilities. The modular platform allows for further integration into an automated work cell for throughput growth or walk-away time.
Imaging	White light and fluorescence	White light and fluorescence	White light and fluorescence	White light and fluorescence
Picking capacity	3000 colonies per hour in white light, 2000 colonies per hour in fluorescent light	3000 colonies per hour in white light, 2000 colonies per hour in fluorescent light	3000 colonies per hour in white light, 2000 colonies per hour in fluorescent light	1500 colonies per hr white light and 1000 colonies per hour in fluorescence mode
Colony selection criteria	Size, proximity, roundness, fluorescence intensity	Size, proximity, roundness, fluorescence intensity	Size, proximity, roundness, fluorescence intensity	Size, proximity, roundness, fluorescence intensity
Picking and regional picking	•	•	•	•
Barcode tracking	•	•	•	•
Re-arraying and replication	•	•	•	•
Gridding	•	•	•	•
Plating and streaking		Only QPix 460	•	
Agar to agar			•	
Robotics integration			•	
Shaking incubator			•	
Liquid handler			•	
PCR			•	
Sealer/peeler			•	•
ELISA			•	•
Destination plate capacity	Picking: 12 plates Replicating and re-arraying: Maximum of 20 plate positions	QPix 450: Up to 156 standard SBS plates, 52 standard SBS plates per stacker, up to 3 stacker lanes. QPix 460: Up to 104 standard SBS plates, 52 standard SBS plates per stacker, up to 2 stacker lanes.	Configurable and expandable with automation. No limit to number of plates.	Picking: 4 plates Replicating and re-arraying: Max – 8 plates
Stackers		2 or 3 Stacker Lanes	Plate Hotels	
Source plate capacity	Without manual intervention: 1 x 15 cm petri dish; 5 x 9 cm petri dishes; 2 x OmniTrays; 1 x 22 cm QTrays	Without manual intervention: 2 x 15 cm petri dish; 10 x 9 cm petri dishes; 4 x OmniTrays; 2 x 22 cm QTrays	Automation mode: 1-well Omnitrays, 8-well Omnitrays Manual mode: Qtrays, Petri Dishes, Omnitrays, SBS plates	Manual Mode: 1 x 15 cm petri dish; 5 x 9 cm petri dishes; 2 x OmniTrays; 1 x 22 cm QTrays Automation mode: 1-well Omnitrays, 8-well Omnitrays
Walkaway time	25 minutes at a time – only return to swap destination plates after 12 are full	QPix 450: 156 plates x 96 colonies per plate = 14,976 colonies picked in 4.5 hours QPix 460: 104 plates x 96 colonies per plate = 9,984 colonies picked in 3 hours	Entire duration of run	



대표전화 02-3471-9531
 학술문의 엠디서비스.com
 주소 서울특별시 강남구 테헤란로 623 삼성빌딩 15층