

Your complete solution for sample prep in process development and protein characterization.

The JANUS™ G3 BioTx™ automated workstations enable consistent small-scale protein purification and sample prep for analytical protein characterization required to support quality by design experimentation in both upstream and downstream processes. Automating sample prep provides researchers more time to focus on new analytical tests obtaining critical information earlier in the protein development pipeline. Time and labor savings accelerate project workflows and thus, commercialization of protein therapies.

Automate four modes of small-scale protein purification on a single platform

- PhyNexus® PhyTips® columns
- GE® PreDicator™ plates/ Pall® AcroPrep™ filter plates
- OPUS® RoboColumns® Chromatography Columns
- GE® His GraviTrap™ columns

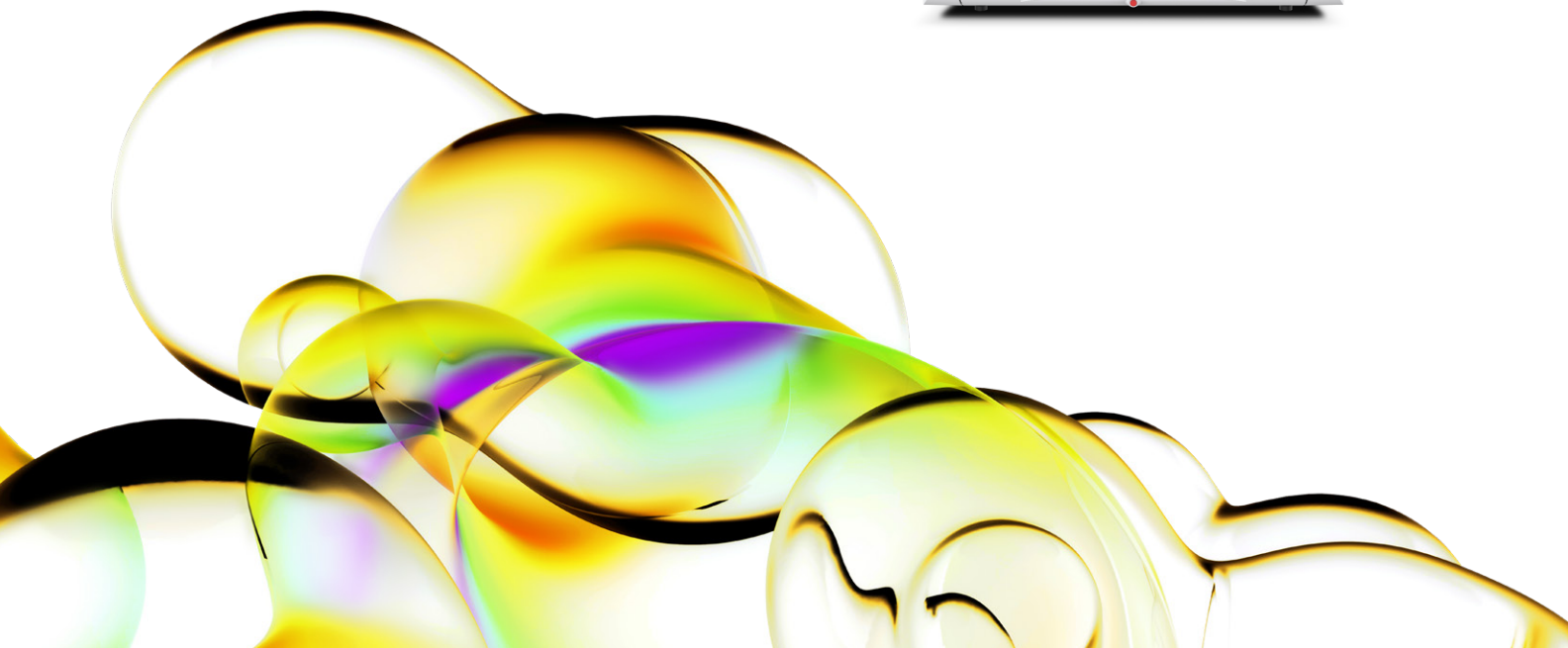
Consistent, time-savings results for your biotherapeutics development:

- Expression optimization
- Cell line selection
- Purification process development
- Bioanalytical sample prep

JANUS™ G3 BioTx™ Workstations



For research use only. Not for use in diagnostic procedures.



Maximize productivity with automated purification methods

The JANUS G3 BioTx workstations accommodate column, tip, and batch chromatography modes on one platform. With just one instrument, you can achieve rapid analysis of

a range of sample volumes and concentrations to support Quality by Design experimentation in both upstream or downstream processes.

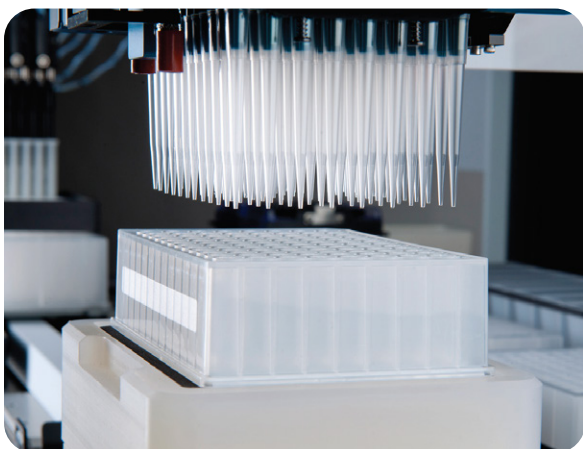
A



PhyNexus® PhyTips® Columns

- Purify up to 96 samples in 15 min
- Column bed sizes from 5-160 μ L volumes
- Achieve high sample binding capacity and purity by passing samples back and forth through column
- Highly purified products with high concentrations achieved with low elution:bed volume ratio

B



GE® Predictor® Plates/Pall® AcroPrep™ Filter Plates

- Rapidly screen binding, wash and elution conditions in 96 well plate formats
- Use centrifugation or vacuum filtration (optional gripper for full automation)
- Explore design space to optimize process development using DoE

C

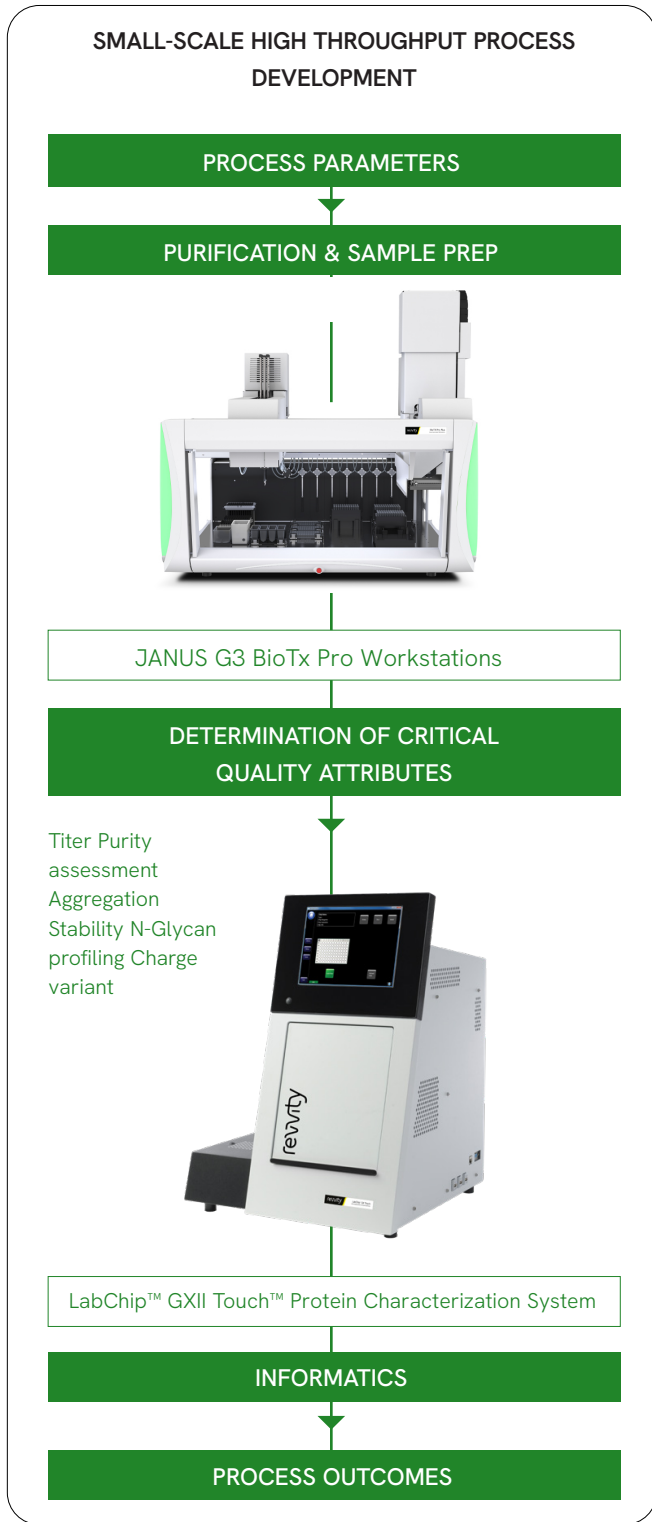


OPUS® RoboColumns® Chromatography Columns

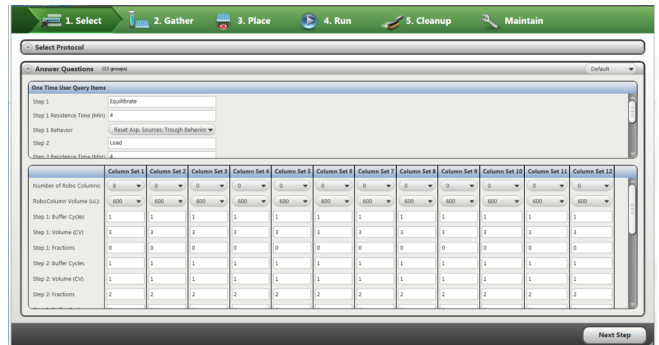
- Miniaturized column chromatography predictive of process scale-up
- Purify 96 samples/day or more depending upon load volume
- 50, 100, 200 and 600 μ L bed sizes; variety of resin types
- Sample volumes from 100 μ L up to 48 mL or more
- Automate sample loading and fraction collection; adjustable height accommodates multiple collection plates

Figure 1. The JANUS G3 BioTx workstation accommodates (A) PhyNexus® PhyTips® columns, (B) filter plates and (C) OPUS® RoboColumns® Chromatography Columns on one platform. Purification methods include affinity, ion exchange, reverse phase and gel filtration.

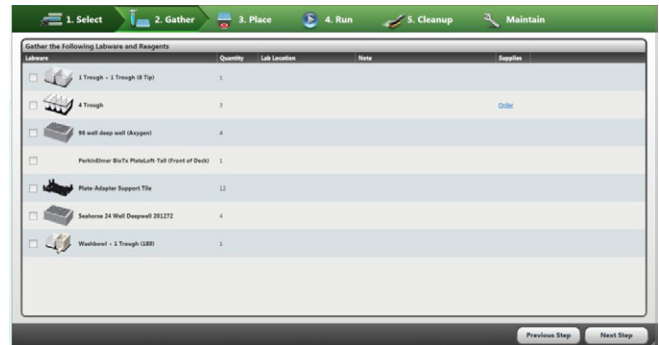
Perform Qbd analysis in hours, not weeks



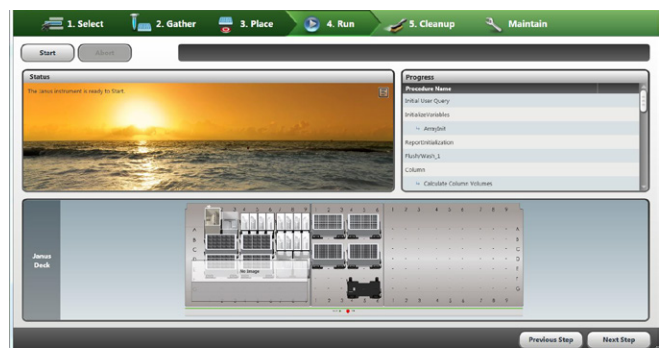
Simple to use protocols—designed for the scientist



Quickly enter experiment variables to easily translate your experimental design



Easily confirm placement of samples and labware on the workstation deck (top) and view workstation status or setup email confirmation for when process is complete (bottom).



Wizard driven JANUS G3 application assistant guides preparation and execution of small-scale purification and sample prep

- Enables true walk-up usability without need for advanced training
- Flexibility to easily modify purification parameters

Figure 2. With the JANUS G3 BioTx™ workstations and LabChip GXII Touch protein characterization system, you can purify and analyze in one day what typically takes weeks with alternative methods. Explore a broader range of experimental conditions and save in both the development time and cost of your biopharmaceuticals research.

Choose the JANUS G3 BioTx Pro or JANUS G3 BioTx Pro plus workstation depending upon your throughput and capacity needs

- Eight tip independent pipetting arm designed for multiple protein chromatography applications:
PhyNexus® PhyTips® columns, GE® Predictor® Plates/
Pall® AcroPrep™ filter plates, OPUS® RoboColumns®
chromatography columns, GE® His GraviTrap™ columns

JANUS G3 BioTx pro workstation

- 12 deck positions
- Compact footprint



- Flexible control software and wizard driven application assistant
- Automated fraction collection station with housing for up to 96 small scale columns

JANUS G3 BioTx pro plus workstation

- 24 deck locations
- MDT 96-tip head for parallel purification methods
- Integrated gripper for labware positioning
- Integrated vacuum manifold



Complete your small-scale protein purification workflow with the LabChip GXII Touch protein characterization system

The LabChip GXII Touch protein characterization system complements QbD and small-scale protein purification studies by offering high throughput, consistent and precise analysis of protein samples.

- Characterize seven protein critical quality attributes with comparable performance to traditional CE
- At 70X greater throughput
- And 75% greater labor savings



For more information, visit www.revvity.com

revvity