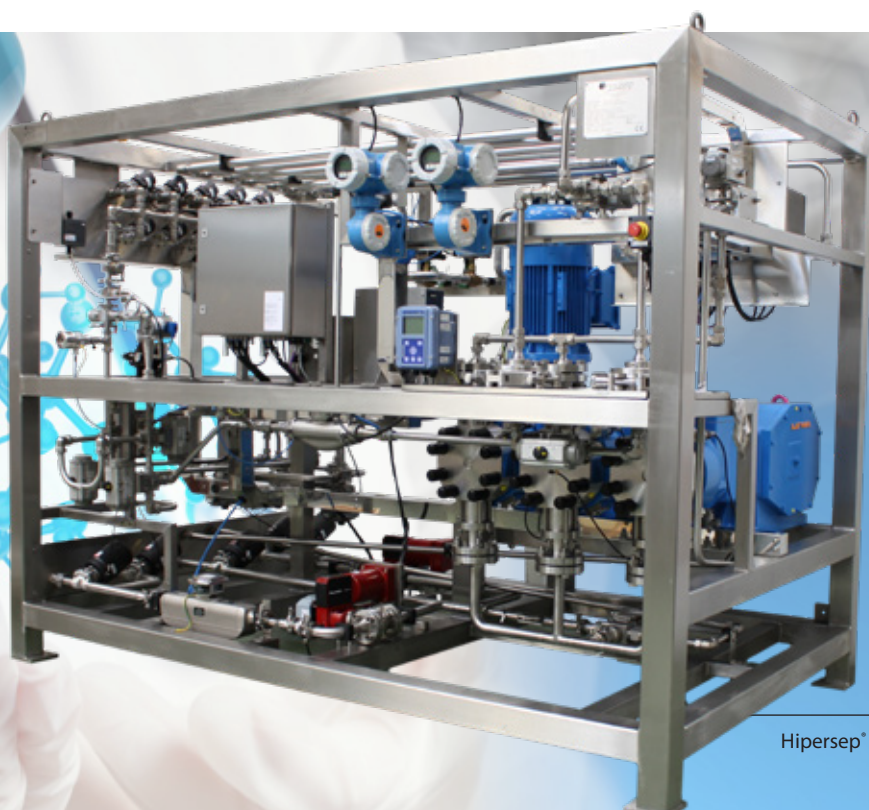


HIPERSEP[®] skids

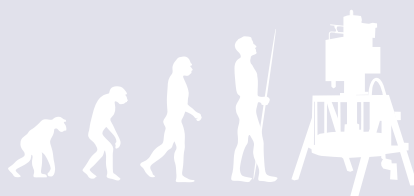
PREPARATIVE AND INDUSTRIAL HPLC SOLUTIONS



Hipersep[®] XXL



- High performance, high reproducibility
- Accurate and consistent gradient
- Automated automation for linear and easy scale-up
- Explosion proof design (ATEX & Class Div)
- GAMP, FDA 21CFR Part 11 validated, simple and intuitive software
- Full regulatory compliance (FAT/SAT, IQ/OQ)



Best-in-Class HPLC skids for pharmaceutical applications

With a strong track record in developing, using and supplying preparative HPLC solutions, Novasep offers state-of-the-art Hipersep® skids for intermediate purification and polishing of APIs, biologics and other small molecules (peptide, oligo, insulin...), from development to large-scale commercial phase.

Get the best from chromatography with Hipersep® skids



Control

Ensure high reproducibility with our smart control system.

Sanitary

Easy-to-clean, minimized risk of cross-contamination.

Saving

Save time and solvent consumption with our accurate gradient module (<1%).

Compact

Fits anywhere!



Flexible

Over a 1,000 possible configurations fit with columns from 50 to 150mm I.D.

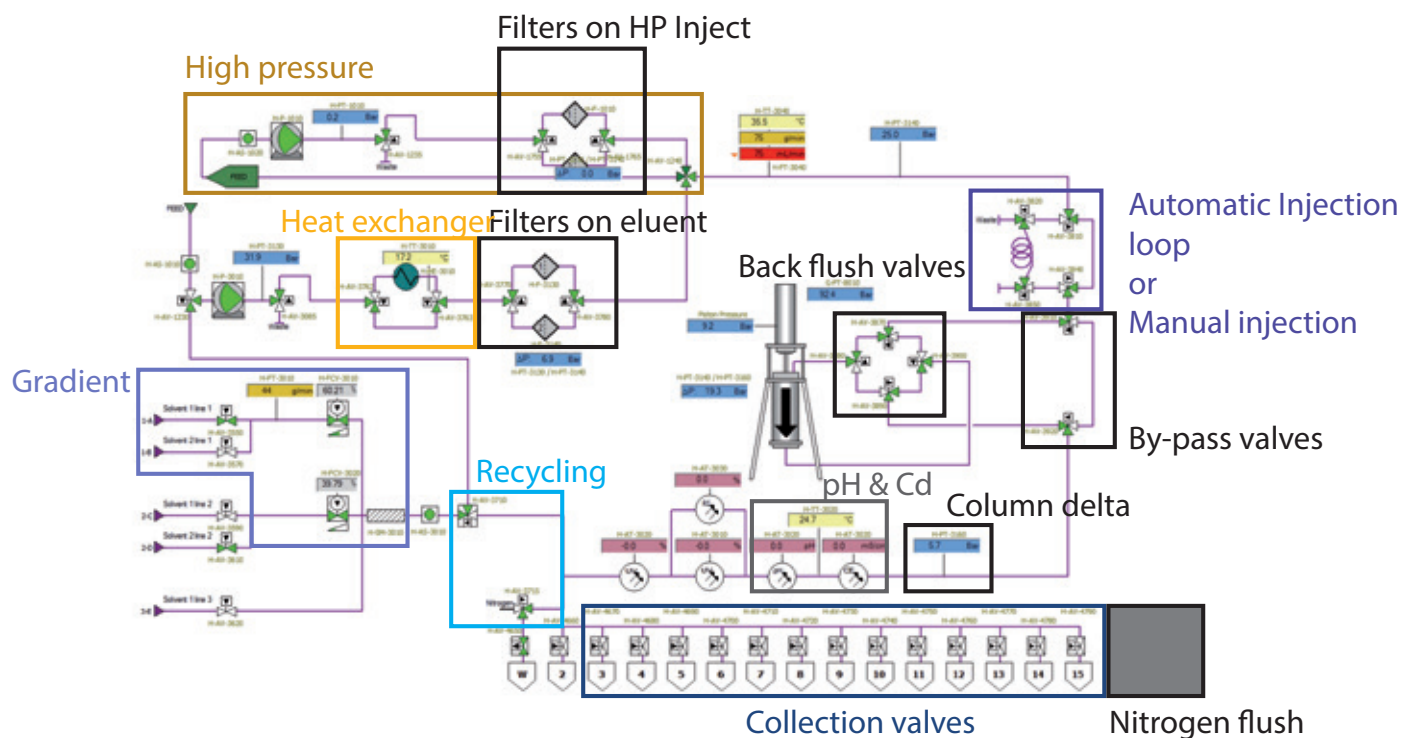
Cyclojet

Increase productivity of poor resolution processes.

Design

Save on your skid maintenance costs through our robust and industrial design made for 24/7 operations.

Design your own skid with our modular concept



Whatever your scale, Novasep can provide the right solution

From lab to manufacturing scale operations

Hipersep®	Pilot	Process M	L	XL	XXL	Above
Productivity range (kg/day) in Batch mode	0,2-4	2-33	17-57	29-106	53-165	
Productivity range (kg/day) in Cyclojet	8-12	14-99	114-171	212-318	330-495	
Typical Prochrom® column diameter (mm)	50/80/110/150	110/150/200/300/450	450/590	590/800	800/1000	
Skid flowrate (L/h)	6*-90	60-500 / 20-200 (Low Flow version)	300-1000	500-1600	1000-2500	
Gradient accuracy (%)	< 1					
Max pressure (bar)	100				70	
Explosion proof	Atex Zone II Class 1 Division 2					
Certifications / Equipment standards	CODAP/ASME for columns CE/UL for systems					
Configurable options	High pressure injection module, Eluent & feed filters, Heat exchanger, Backflush configuration valves, pH conductivity, Nitrogen flush for product recovery, Up to 10 collections, Solvent recycling, Cyclojet process					

*Gradient accuracy <2% below 10L/h

Simple and intuitive Hipersep® SC Software

21CFR Part 11 & GAMP 5 Ready Control Software
Hipersep® SC makes your chromatography processes user-friendly, robust and traceable:

- Intuitive recipe building
- Manual/automatic operations
- Real-time monitoring of all process parameters
- Several modes of smart, event-based fraction collection
- Batch management and historical data logging

Fast Global Support



Capitalized on Novasep's know-how by becoming an expert on Novasep technologies with our wide range of training programs (equipment, process, maintenance, software).



Through our process development team, we have the possibility to develop, optimize and industrialize your valuable molecules.



Stay up to date with our software upgrade services package.



Maximize your up time by managing your OPEX with different levels of maintenance contracts.



Based on your needs, you have the best of Novasep readily on your site. Hot line available 24 hours a day, 7 days a week, to respond to any urgency.



Services and technologies
for the life science industries

www.novasep.com

Contact our experts at
novasep@novasep.com

EUROPE

Novasep Equipment Solutions
Site Eiffel
81, boulevard de la Moselle - BP 50
54340 Pompey - FRANCE
Phone +33 383 497 000

AMERICAS

Novasep LLC
23 Creek Circle
Boothwyn, PA
19061 - USA
Phone +1 484 361 6000

ASIA

Novasep Asia
Building No.4,
Lane 1690, Zhangheng Road,
Zhangjiang Hi-Tech Park,
Shanghai 201203 - CHINA
Phone: +86 (0)21 6045 1600

INDIA

Novasep India
Four Square Elegant,
Block A, Flat 110,
Madhavapuri Hills, Road 5,
Hyderabad 50050 - INDIA
Phone: +91 9866573377