

Clean In Place System (CIP) – Essential Range

To Be Connected to a Manual Filtration Skid – Essential Range

Product Information

All Manual filtration skids from the Essential range can be connected to a CIP system. It produces and maintains hot water up to 90 °C | 194 °F. It supplies filtered water for the filling line (filtration skid, piping and bottling machine) either for cleaning purposes at 50 °C | 122 °F or for sterilization at 90 °C | 194 °F. Exists in 1000, 1500 and 2000 liters.

Applications

It supplies filtered water for the filling line (filtration skid, piping and bottling machine) either for cleaning purposes at 50 °C | 122 °F or for sterilization at 90 °C | 194 °F. The CIP System optimizes water consumption for the rinsing, cleaning and sterilization phases.



CIP System configuration dependent on filling line specification

CIP System	Up to 2000 Btls/hr		Up to 5000 Btls/hr	
	2 stage filtration	3 stage filtration	2 stage filtration	3 stage filtration
1000 L	•	•		
1500 L			•	
2000 L			•	•

Basic Configuration

- Cleaning flow control by manual valve and hot water pump
- Hot water unit (90 °C | 194 °F) by electric resistor
- Hot water unit (50 °C | 122 °F) by thermostatic mixing valve
- Prefiltration of incoming water
- Water temperature maintained
- Programmable tank filling and water heating to align with off-peak hours

Advanced configuration – available options

- Automated sterilization module with integrated data traceability
- Recirculation loop accessory for filtration skids & bottling lines, limiting water loss from sterilization cycles (up to 80% of recovery)
- Integrity testing automation module for wine final filter on the filtration skid with data traceability
- Automated dosing kit for oenological adjuvants on the filtration skid
- All modules are controlled through PLC (Programmable Logic Controller) System.

Equipment on the skid

	1000 L CIP System	1500 L CIP System	2000 L CIP System
Electrical pump: 10 m3/hour	x	x	x
Electrical frame	x	x	x
PLC option dependent	x	x	
Temperature sensors	x	x	x
Level sensors	x	x	x
Manual and automatic valves. Butterfly and sphere	x	x	x
Tank capacity	1000 L	1500 L	2000 L
Electrical resistance	2 × 12 KW	2 × 18 KW	2 × 18 KW
Water filter	1 × 30"	1 × 30"	1 × 30"

Options

	1000 L CIP System	1500 L CIP System	2000 L CIP System
Automatic start hot water pump	x	x	x
Automatic sterilization	x	x	x
Recirculation	x	x	x
Dosing arabic gum	x	x	x
Integrity test	x	x	x
Recirculation loop	x	x	x

Technical Data

Main construction material	304L stainless steel
Piping surface finish	<ul style="list-style-type: none"> ▪ Inner surface Ra < 0.8 µm / welded seam area Ra < 1.6 µm ▪ Outer surface Ra < 1.0 µm
Insulation specification	50 mm (1.97 in) polyester + external PVC cover
Inlet & outlet connection	DN25
Power supply	Control cabinet: 400 V 3P+PE 40 Kw
Frame design	Square pipes structure
Frame feet	Anti-slip
Frame surface finish	Mechanical polishing

Dimensions and Weight (approx.)

	W × D × H	Weight
1000 L	1250 × 1250 × 2217 mm (49 ³ / ₁₆ × 49 ³ / ₁₆ × 87 ¹ / ₄ in)	630 kg (1,389 lbs)
1500 L	1500 × 1400 × 2221 mm (59 ¹ / ₁₆ × 55 ¹ / ₁₆ × 87 ⁷ / ₁₆ in)	810 kg (1,786 lbs)
2000 L	1500 × 1400 × 2755 mm (59 ¹ / ₁₆ × 55 ¹ / ₁₆ × 108 ⁷ / ₁₆ in)	930 kg (2,050 lbs)

Ordering Information

Product code	Designation
Clean In Place system Essential range	
CIP-ESS-1000	Manual clean in place system 1000L
CIP-ESS-1500	Manual clean in place system 1500L
CIP-ESS-2000	Manual clean in place system 2000L


Clean In Place system Essential range – List of options	
FSKID-ESS-CIP-OPT1	Automated pump start
FSKID-ESS-CIP-OPT2	Automated sterilization program
FSKID-ESS-CIP-OPT3	Automated sterilization program + water recovery
FSKID-ESS-CIP-OPT4	Automated sterilization program + automated integrity test
FSKID-ESS-CIP-OPT5	Automated sterilization program + oenologic product dosage

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com