

CRYOLYS® EVOLUTION



Cooling unit for sensitive sample preparation

- **Efficient cooling:** before, during and after processing
- **No need for specific air pressure or power supply**
- **Accurate temperature** monitored by Precellys® Evolution software
- **Invisible footprint on the bench**



CRYOLYS® EVOLUTION

Keep your thermo-sensitive molecules "in shape"!

Monitored low temperature ensures integrity for specifically targeted molecules like proteins and RNA, giving downstream applications better efficiency/sensitivity.

Cryolys® Evolution is the only patented integrated cooling unit preventing the increase of temperature during lysing. Based on dry ice sublimation, it enhances the efficiency of molecular extraction, resulting in premium quality analysis, maintaining the desired temperature between 0 and 10°C before and during homogenization.

Thanks to its ingenious design, the Cryolys® Evolution is now a real extension of the Precellys® Evolution homogenizer. Positioned on the unit's lid, the system is powered and managed directly by the Precellys® Evolution software, offering an invisible footprint in the lab and a complete control of the temperature during samples grinding.

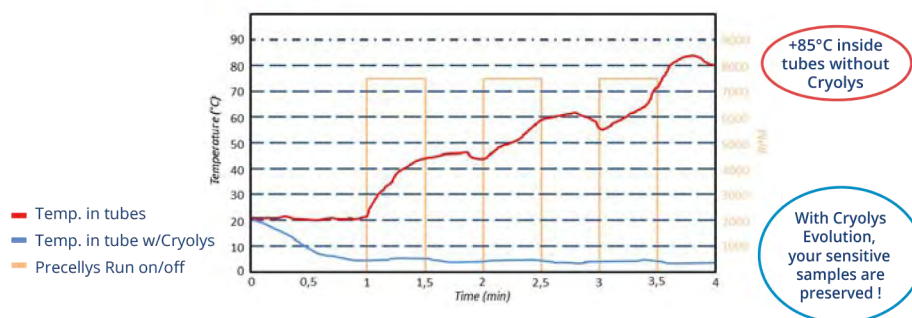
- Achieve higher yield, greater quality and increased functionality
- Protect from enzymatic activities
- Ensure native protein conformation conservation

Cryolys® Evolution - Principle



► Tubes temperature profile with Cryolys® Evolution

7500rpm, 3x30s, 30s breaks



Technical specifications

SIZE:

L: 335mm, W: 335mm, H: 100mm

WEIGHT:

Full unit: 2,6 kg
Empty bowl: 1,3 kg

BOWL CAPACITY: 1,5 kg of dry ice

DRY ICE CONSUMPTION:

1,5 kg for 30 min (run time)

NORMS: CE

NOISE: < 60 dB

PATENTED TECHNOLOGY

REQUIREMENT

Unit plugged to a Precellys® Evolution

CATALOG NUMBER

EQ05648-300-RD001.0

Precellys homogenizers range

Discover the full range of products



Precellys® Evolution
Super Homogenizer



Minilys
for personal use



Precellys® 24
for high throughput



Lysing
kits consumables