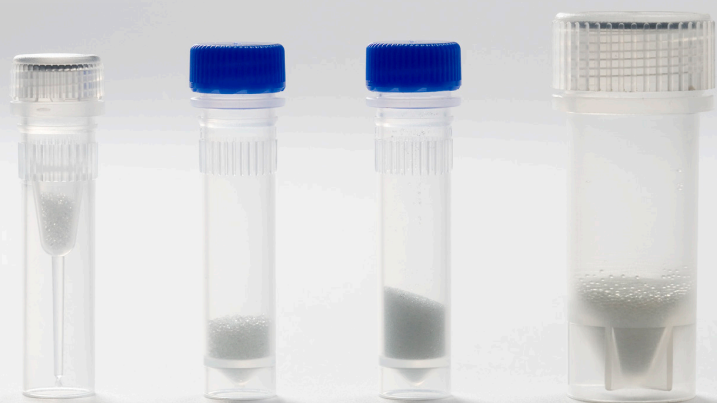


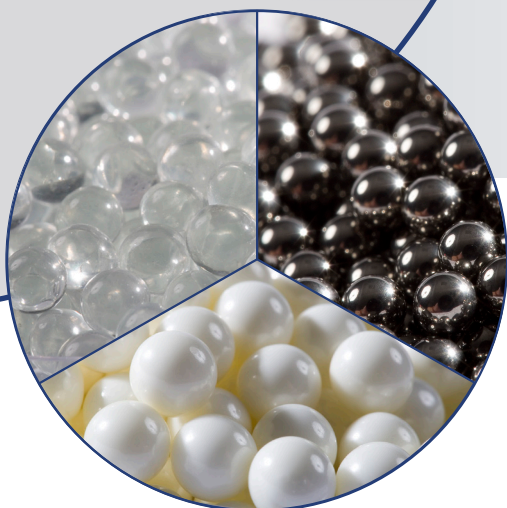


Precellys

The Science of Lysing



0,5ml tubes
2ml tubes
7ml tubes



Precellys Lysing kits for tissue homogenizer

Precellys Lysing Kits are prefilled tubes with specific beads and are suitable for use with all biological samples prepared for DNA, RNA, protein or drug extraction.

- More than 15 references for soft samples to hard samples
- 0,5ml – 2ml – 7ml ready-to-use tubes
- Wide range and mix of beads for optimized results
- Full range of ceramic beads

1000 protocols
available on
www.precellys.com

www.precellys.com

bertin
TECHNOLOGIES

A company of **CNIM** Group

PRECELLYS LYSING KITS

A complete range for biological samples preparation

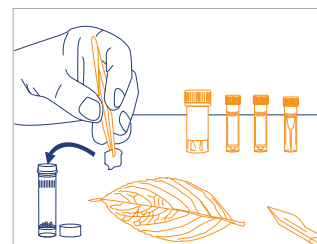


The Science of Lysing

Precellys LYSING KITS

Enjoy all the benefits of **Precellys' powerful bead-beating technology** with the 15 available references of Precellys Lysing Kits! Precellys Lysing Kits are robust tubes prefilled with high quality beads and manufactured under clean-room environment. **Optimize your preparation of biological samples:** animal tissue, human tissue, plant tissue, microorganisms... whatever the precious sample you wish to prepare, find a suitable Precellys Lysing kit from the **15 available references:** choose the right **tube size** (0.5ml, 2ml or 7ml), **bead type / diameter** (ceramic, glass, stainless steel ...) and you will be sure to obtain **excellent DNA, RNA, Protein and Drug recovery!**

- Save time and improve your lab practices by following tried & tested scientific publications: use our web-based Application Center to find the right Lysing Kit for your sample: just enter your sample name and get access to examples, lysing kits and protocols that current Precellys' users are already using!



Animal and Human tissue homogenization

Brain, liver, heart, muscle, skin, kidney, lung, spleen, colon, faeces etc. Also pulverization of difficult tissue samples such as bones, hair, cartilage, etc.

Whole organs can even be prepared thanks to 7ml Lysing Kits.



Plant tissue grinding

Leaves, oats, tomato seeds, cotton, soil etc.

Also pulverization of hard materials such as seeds, roots, rice etc.

Large samples can also be prepared thanks to 7ml Lysing Kits.



Micro-organism lysing

Small beads facilitate the crushing of micro-organism walls.

Yeast, fungi, spores, mould, bact. gram+/gram-, mycobacteria, cyanobacteria, algae etc.

0.5ml Lysing Kits are especially suited to small sample amounts.



- Log on now to www.precellys.com to find the most suitable lysing kit for your sample!

Precellys range

- Precellys Lysing Kits are specially designed for use with the entire Precellys range:



Precellys 24
for high throughput



Precellys 24-Dual
for flexibility



Cryolys
for sensitive sample



Minilys
for personal use

Technical specifications

QUANTITY OF TUBES PER KIT: 50

TUBES & CAPS MATERIALS:
Polypropylen Co-Polymered

TEMPERATURE RANGE OF USE:
-80°C to 100°C

AUTOCLAVE: standard procedure

CHEMICAL COMPATIBILITY: compatible
with most of extraction buffer (Acid
Phenol, Guanidine thiocyanate,
acetonitril etc.)

LOADING CAPACITY 0.5ml: from 1mg
to 30mg tissue or 5µL to 200µL culture

LOADING CAPACITY 2ml: from 30mg
to 1g tissue or 200µL to 1.6ml culture

LOADING CAPACITY 7ml: from 30mg
to 5g tissue or 200µL to 6ml culture

COMPATIBLE WITH PRECELLYS 24,
PRECELLYS 24-DUAL, CRYOLYS
AND MINILYS.

1000 protocols
available on
www.precellys.com

Find your distributor on www.precellys.com

BERTIN TECHNOLOGIES

FRANCE

Phone: +33 139 306 169

BERTIN Corp.

MD, USA

Phone: +1 240 428 1047

precellys@bertin.fr