

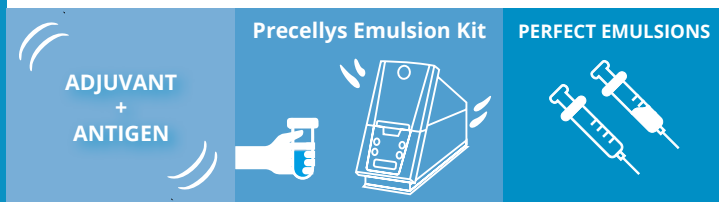
Precellys Emulsion Kit

Get perfect emulsions



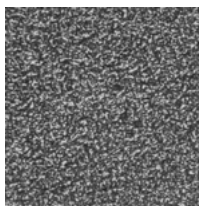
- **Easy:** put your Adjuvant and Antigen (peptide, bacteria...) solutions in the tube and let the instrument mix for you
- **Fast and Efficient:** only 2 min shaking time on Minilys or Precellys Evolution for a perfect emulsion
- **Flexible:** prepare 2 to 8mL of emulsion
- **Safe & Easy:** no risk of accidental exposure thanks to the single tube concept and no Repetitive Stress Injuries (RSI)

An efficient method...

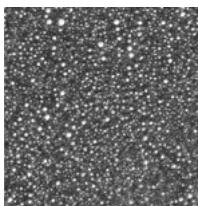


...to prepare homogenous and stable emulsions

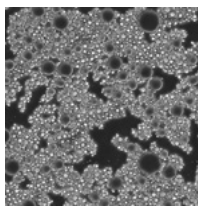
Precellys Emulsion kit



2-syringe method



Vortex



Precellys Emulsion Kit can be used for:

- Preparation of emulsions with Freund's Adjuvant + Antigens
- Monoclonal & polyclonal antibody production
- Preclinical animal models for research & in industry
- DTH experiments (Delayed Time Hypersensitivity)

REFERENCE: D34200

Content

- 10 autoclaved pouches including:
- 1x Emulsion tube and cap
 - 1x specific Emulsion tube plunger

DISCOVER OUR LIFE SCIENCES PRODUCT RANGE



AIR SAMPLING



SAMPLE
PREPARATION



ASSAY KITS



CELL IMAGING
SYSTEM

Bertin Technologies - Head Office

10 bis Avenue Ampère - Parc d'activités du Pas du Lac
78180 Montigny-le-Bretonneux FRANCE