

PRODUCT NAME

15 ml Bioruptor® Tubes

Cat. No: C30010017**Format:** 50 rxns**Description**

The 15 ml Bioruptor® Tubes strongly improve protein extraction efficiency. The tubes have been thoroughly validated for use on the Bioruptor® Pico (B01060001) and are strongly recommended for extraction of proteins using this instrument.

Characteristics

- Clear
- Steril