

PRODUCT NAME

Protein Extraction Beads

Cat. No: **C20000021**

Format: 20 g

Description

The protein extraction beads have been extensively validated to be used in combination with Diagenode's Bioruptor® Standard, Plus and Pico to ensure efficient disruption and homogenization of fresh and frozen tissues. The Protein Extraction Beads guarantee high yields of proteins.

Storage & Stability

Store at room temperature.

Precautions

This product is for research use only. Not for use in diagnostic or therapeutic procedures.