

CUTANA™ CUT&RUN 8-strip 0.2 mL tubes

| | | | |
|-------------------|-------------|------------------|------------|
| Catalog No | 10-0009 | Pack Size | 120 strips |
| Lot No | 23146005-02 | | |

DESCRIPTION

CUTANA™ CUT&RUN 8-strip 0.2 mL tubes and attached flat caps are made from high quality resins without lubricants or additives for reproducible results. The tubes are autoclavable and supplied RNase, DNase, and DNA free.

FEATURES

These clear tubes are supplied in boxes of 120 strips, compatible for 960 CUTANA™ CUT&RUN reactions.

APPLICATION NOTES

Suitable for cell-based sorting and molecular biology applications, including CUTANA™ CUT&RUN and CUT&Tag applications. See the CUTANA™ assay protocols for more details (epicypher.com/protocols).

PRODUCT IMAGES



FIGURE 1 CUTANA™ CUT&RUN 8-strip 0.2mL tubes. 8-strip 0.2 mL tubes enable multichannel pipetting for improved sample handling and increased experimental throughput in CUT&RUN applications. The tubes are pictured with CUTANA™ Concanavalin A (ConA)-Conjugated Paramagnetic Beads (EpiCypher 21-1401), which are used for immobilizing cells or nuclei during the CUT&RUN procedure. CUTANA™ Concanavalin A-Conjugated Paramagnetic Beads are not included in the purchase of this product.



FIGURE 2 CUTANA™ CUT&RUN 8-strip 0.2mL tubes. 8-strip 0.2 mL tubes are compatible with the Magnetic Separation Rack (EpiCypher 10-0008) for use during all bead separation steps in the CUT&RUN procedure, including sample immobilization with ConA beads and library preparation. Magnetic Separation Rack is not included in the purchase of this product.