

CUTANA™ 8-strip 0.2 mL tubes

Catalog No	10-0009	Pack Size	120 strips
Lot No	25062009-01		

DESCRIPTION

CUTANA™ 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 or CUT&Tag 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™ 8-strip 0.2 mL tubes. 8-strip 0.2 mL tubes enable multichannel pipetting for improved sample handling and increased experimental throughput in CUT&RUN and CUT&Tag 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 and CUT&Tag procedures. CUTANA™ Concanavalin A-Conjugated Paramagnetic Beads are not included in the purchase of this product.



FIGURE 2 CUTANA™ 8-strip 0.2 mL 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 and CUT&Tag procedures, including sample immobilization with ConA beads and library preparation. Magnetic Separation Rack is not included in the purchase of this product.