

Magnetic Separation Rack, 0.2 mL Tubes

| | | | |
|-------------------|-------------|------------------|--------|
| Catalog No | 10-0008 | Pack Size | 1 Rack |
| Lot No | 25055017-81 | | |

DESCRIPTION

Magnetic Separation Rack for use with up to sixteen 0.2 mL PCR tubes for separation of paramagnetic beads in a variety of cell-based or molecular biology applications, including Cleavage Under Targets and Release Using Nuclease (CUT&RUN) and Cleavage Under Targets and Tagmentation (CUT&Tag).

FEATURES

Solid aluminum construction with hard coat anodizing for durability.

Strong permanent Neodymium magnets for fast and effective bead separation.

Compatible with standard 0.2 mL PCR tubes, either as individual tubes or strip tubes.

Double-sided for simultaneous processing of up to 16 samples.

Can be used on either small volumes (5 - 50 μ L, Figure 1) or larger volumes (50 - 200 μ L, Figure 2).

APPLICATION NOTES

Suitable for cell-based sorting and molecular biology applications using paramagnetic beads, including CUTANA™ CUT&RUN and CUT&Tag applications. Also needed for the CUT&RUN Library Prep Kit (epicypher.com/14-1001).

For more information, see our CUTANA protocols (epicypher.com/protocols).

PRODUCT IMAGES



FIGURE 1 Magnetic Separation Rack, front view. Note that the 5 - 50 µL is up, and that the magnets are toward the bottom of the rack, so that the beads will be sorted near the bottom of the tubes.



FIGURE 2 Magnetic Separation Rack, oblique view. Rack shown with PCR strip tubes and paramagnetic beads. Note that the 50 - 200 µL configuration is being used, and that the magnets are sorting the beads to the side of the tubes.