

## Magnetic Separation Rack, 1.5 mL Tubes

|                   |             |                  |        |
|-------------------|-------------|------------------|--------|
| <b>Catalog No</b> | 10-0012     | <b>Pack Size</b> | 1 Rack |
| <b>Lot No</b>     | 25266011-81 |                  |        |

### DESCRIPTION

Magnetic Separation Rack for use with up to six 1.5 mL microfuge 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 1.5 mL microfuge tubes.

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

### APPLICATION NOTES

Suitable for cell-based sorting and molecular biology applications using paramagnetic beads, including CUTANA™ CUT&RUN and CUT&Tag applications. For more information, see our CUTANA protocols ([epicypher.com/protocols](http://epicypher.com/protocols)).

### PRODUCT IMAGES



**FIGURE 1 Magnetic Separation Rack, front view.** The rack is double-sided and can hold up to six 1.5 mL microfuge tubes at a time.