

# Magnetic Separation Rack, 1.5 mL Tubes

**Catalog No.** 10-0012

**Lot No.** 21068003-02

**Unit Size** 1 Rack



**EpiCypher**®

## Product 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, with two sets of magnets.

## Application Notes:

Suitable for cell-based sorting and molecular biology applications using paramagnetic beads, including CUTANA ChIC/CUT&RUN and CUT&Tag applications. Also useful for NGS library preparation.

See detailed CUTANA assay protocols here:

<https://www.epicypher.com/resources/protocols/>

## References:



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

This product is for *in vitro* research use only and is not intended for use in humans or animals.

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • [www.epicypher.com](http://www.epicypher.com)