

## MagniSort™ Magnet

Catalog Number: MAG-4902

**RUO: For Research Use Only. Not for use in diagnostic procedures.**

### Product Information



**Contents:** MagniSort™ Magnet

**Catalog Number:** MAG-4902

**Handling Conditions:** This magnet is comprised of strong rare earth magnets. Rare earth magnets are sensitive to heat above 175°F (80°C). Keep away from extreme heat source.



**Batch Code:** Refer to vial



**Use By:** Refer to vial



**This magnet generates a strong magnetic field. Keep away from pacemakers, credit cards, magnetic I.D. cards, watches, computer monitors and hard disks to prevent damage to these devices.**

### Description

The MagniSort® Magnet, 5 mL, is designed to be used with 12 x 75 mm round bottom polystyrene tubes, and can be used to magnetically separate up to  $2 \times 10^8$  cells. For detailed instructions, follow the protocols provided with each separation kit.

### Applications Reported

The MagniSort® Magnet, 5 mL, has been reported for use in magnetic cell separation.

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • [thermofisher.com/ebioscience](http://thermofisher.com/ebioscience) • [info@ebioscience.com](mailto:info@ebioscience.com)