

Description

This Cell Isolation Magnetic Tube Rack is designed to accompany our magnetic cell isolation products, such as the CD19 Positive Cell Isolation Kit (BPS Bioscience #78564). This magnet can quickly separate cells labeled with magnetic beads and is designed to allow users a near 360° view of their cell samples during separation. This magnet can accommodate both 15 and 50 ml tubes simultaneously, which allows flexibility for scaling up or down your isolation samples. Additionally, the rack is compatible with both conical and round-bottom tubes. The compact nature of this magnet allows it to fit easily in various incubators or refrigerators to match desired conditions.



Figure 1: Cell Selection Magnet

Background

Isolating specific cell types from a heterogenous populations such as blood, tissues and cell culture can be time-consuming and laborious and places cells under a great amount of stress. For sensitive populations such as PBMCs, a quick and easy isolation method is ideal to preserve cell viability and function. Magnetic cell isolation removes the stress of repeated centrifugations and allows users to perform separations gently and quickly. This magnet is compatible with all BPS Bioscience Magnetic Cell Isolation Kits.

Applications

Positive and Negative Cell Isolation

Storage Conditions

Store magnet at room temperature away from items sensitive to magnetic fields.

Safety

This product is for research purposes only and not for human or therapeutic use.

This product has a strong magnetic field. Keep away from pacemakers, magnets, computer disks, watches, and other objects that are sensitive to magnetic fields.

Care



After use, wipe the magnet clean with distilled water and then with 70% alcohol to dry the magnet. To disinfect, wipe with 70% alcohol. Do not submerge the magnet in water or liquid.

Troubleshooting Guide

For all further questions, please email support@bpsbioscience.com

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
CD19 Positive Cell Isolation Kit	78564	1 x 10 ⁸ or 1 x 10 ⁹ Cells
5x Cell Isolation Buffer	78563	50 ml