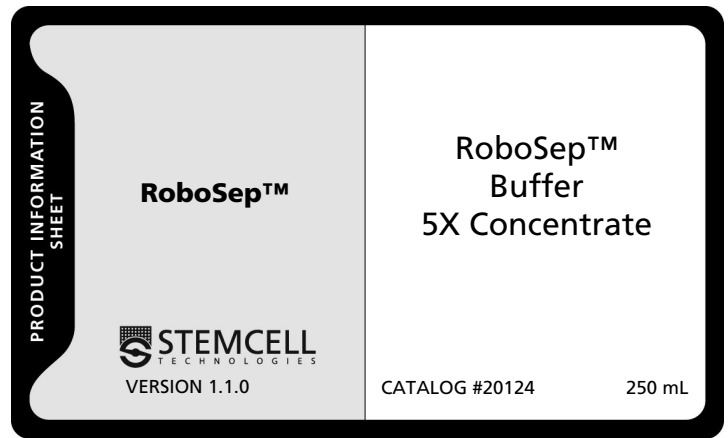


PRODUCT DESCRIPTION

1X diluted buffer from RoboSep™ Buffer 5X Concentrate is recommended for use in EasySep™ cell separation protocols as performed on the RoboSep™ platform. It is also recommended for use in manual EasySep™ cell separation protocols.



COMPONENTS

- 5X Dulbecco's Phosphate Buffered Saline (D-PBS)
- Fetal Bovine Serum (FBS), 10% v/v
- Ethylene Diamine Tetra-acetic Acid (EDTA), 5 mM

The pH range is 7.2 - 7.6.

Product is aseptically manufactured using tightly controlled processes and is sterility tested.

STABILITY AND STORAGE

RoboSep™ Buffer 5X Concentrate: Product stable at 2 - 8°C until expiry date as indicated on label. Avoid repeated freeze-thaw cycles.

1X diluted RoboSep™ Buffer: Product stable at 2 - 8°C for 6 months.

See below for preparation instructions for 1X diluted RoboSep™ Buffer.

DIRECTIONS FOR USE

Prepare 1X diluted RoboSep™ Buffer by adding 1 part RoboSep™ Buffer 5X Concentrate to 4 parts distilled or Type 1 water. Mix well and filter sterilize using a 0.2 micron low protein-binding filter. Dispense into sterile containers.

Refer to the RoboSep™ Technical Manual (Catalog #28940) for directions for use.