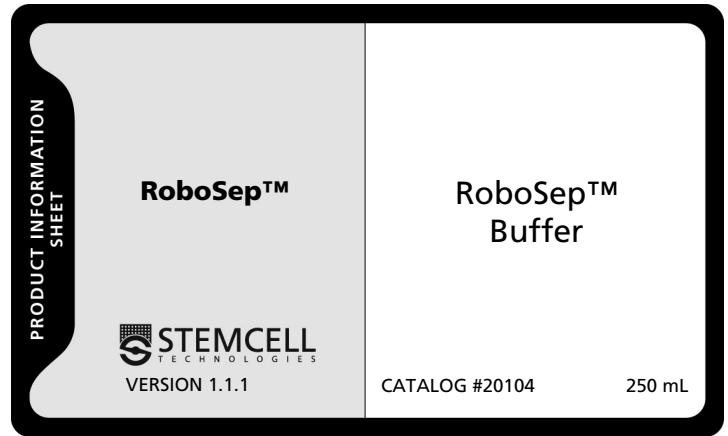


## PRODUCT DESCRIPTION

RoboSep™ Buffer 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

- Dulbecco's Phosphate Buffered Saline (D-PBS)
- Fetal Bovine Serum (FBS), 2% v/v
- Ethylene Diamine Tetra-acetic Acid (EDTA), 1 mM

The pH range is 7.2 - 7.6.

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

## STABILITY AND STORAGE

Product stable at 2 - 8°C until expiry date as indicated on label. Avoid repeated freeze-thaw cycles.

## DIRECTIONS FOR USE

This product has been optimized for use with certain EasySep™ cell separation kits. Please refer to the relevant EasySep™ Product Information Sheets for detailed instructions.

Refer to the RoboSep™ Technical Manual (Catalog #28940) for more information.