

DNASE I SOLUTION

DEOXYRIBONUCLEASE I SOLUTION

Catalog # 07900

1.0 mL/tube

RECOMMENDED FOR:

Addition of DNase I may reduce or prevent clumping of concentrated and/or cryopreserved hematopoietic cell suspensions following thawing.

PRODUCT DESCRIPTION:

This product contains:

Bovine Pancreatic Deoxyribonuclease I
1 mg/mL in Phosphate Buffered Saline

Activity range 2000 - 3000 Kunitz units per mg of protein. One Kunitz unit will produce an A_{260} of 0.001 per minute per mL at pH 5.0 at 25°C, using DNA as a substrate, with a Mg^{++} concentration of 4.2 mM.

STABILITY/STORAGE:

Stable for at least 1 year at -20°C.
Storage at 4°C is not recommended.

Contents are guaranteed sterile if seal has not been tampered with.

DIRECTIONS FOR USE:

1. Thaw cells quickly in a 37°C water bath and transfer cells into a 15 mL or 50 mL sterile tube containing 0.25 to 0.5 mL of 1.0 mg/mL DNase I.
2. Slowly add 10-15 mL of Phosphate Buffered Saline (PBS) or Hank's Balanced Salt Solution (without Ca^{++} or Mg^{++}) dropwise while gently swirling the tube.
3. Fill tube with PBS or Hank's (without Ca^{++} or Mg^{++}) and mix with sterile pipette or by inversion. Do not vortex.
4. Centrifuge at 300 g for 6-10 minutes.
5. After initial cell wash, resuspend cells in 0.1 mg/mL of DNase I. For StemSep™ cell separation procedures, cells can be resuspended to desired cell concentration with the addition of 0.1 mg/mL of DNase to prevent cell clumping during separation. For hematopoietic colony assays, wash cells twice before inoculation into methycellulose-based medium.

**THIS REAGENT IS FOR LABORATORY USE ONLY.
IT IS NOT TO BE ADMINISTERED TO HUMANS.**