

PRODUCT DESCRIPTION

RosetteSep[®] DM-L is a density separation medium designed specifically for use with RosetteSep[®] cocktails for the enrichment of specific human lymphocyte populations (CD3⁺ T cells, CD4⁺ T cells, CD8⁺ T cells, and B cells) from whole blood. The number of desired cells recovered is up to 30% higher when using RosetteSep[®] DM-L instead of Ficoll-Paque[™] PLUS (Catalog #07907/07917/07957/07967).

FORMULATION

- Iodixanol
- Hetastarch
- Sodium Chloride
- Sodium Lactate
- Dextrose
- Calcium Chloride
- Potassium Chloride
- Magnesium Chloride
- Water

Density: 1.081 g/mL

Osmolarity: 290 mOsm

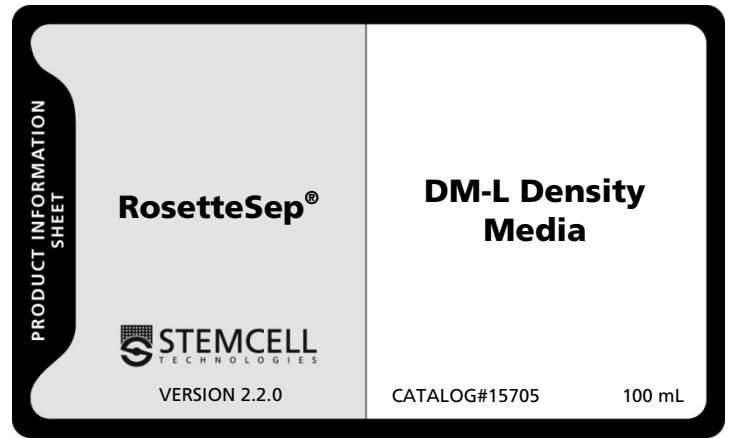
STABILITY AND STORAGE

Store at room temperature (15 - 25°C). Storage at 2 - 8°C is acceptable, but ensure that the medium equilibrates to room temperature (15 - 25°C) and invert bottle to mix contents before use. Keep protected from direct light. Product stable at 15 - 25°C until expiry date as indicated on label. Contents have been sterility tested.

RELATED PRODUCTS

PRODUCT	CATALOG #
<i>Included in kit for:</i>	
RosetteSep [®] HLA Lymphoid Cell Enrichment Kit	15271HLA
<i>Recommended for use with:</i>	
RosetteSep [®] Human T Cell Enrichment Kit	15021/15061
RosetteSep [®] Human CD4+ T Cell Enrichment Kit	15022/15062
RosetteSep [®] Human CD8+ T Cell Enrichment Kit	15023/15063
RosetteSep [®] Human B Cell Enrichment Kit	15024/15064
RosetteSep [®] Human Total Lymphocyte Enrichment Kit	15223/15263

Ficoll-Paque[™] PLUS is a registered trademark of GE Healthcare Ltd.



DIRECTIONS FOR USE

For a list of RosetteSep[®] cell enrichment kits this product can be used with, please see the "Related Products" section on this Product Information Sheet. Please refer to the relevant RosetteSep[®] Product Information Sheets for detailed instructions.

Note: Should not be used with RosetteSep[®] Human Monocyte (Catalog #15028/15068) or RosetteSep[®] Human Progenitor (Catalog #15276, 15026/15066, 15126/15166, 15027/15067 or 15128/15168) Enrichment Cocktails. The density of RosetteSep[®] DM-L is not optimized for the isolation of these cell types.