

PRODUCT INFORMATION SHEET
CD34+ Hematopoietic Stem Cells From Human Umbilical Cord Blood (Frozen)

For Research Use Only

Product Description

- CD34, a transmembrane phosphoglycoprotein, is an established marker of hematopoietic stem and progenitor cells and are used in research areas such as genome editing, immune regulation, hematopoietic reconstitution and regenerative medicine.
- Primary human CD34+ cells are isolated from umbilical Cord Blood using positive immunomagnetic techniques.
- Umbilical cord blood is collected using IRB protocols and/or in accordance with local, state and federal requirements.
- Umbilical cord blood is collected in anticoagulant citrate phosphate dextrose.
- Refer to the lot-specific Certificate of Analyses for product details.

For illustration purposes only. Example order 1×10^6 cells. Actual cell count ordered will be reflected on Certificate of Analysis.

	Specification	Result
Cell Processing Date	Reported	DDMMYYYY
CD34+ Cell Count	$> 1 \times 10^6$ cells	Pass
CD34+ Purity	$> 90\%$	Pass
CD34+ Viability	$> 95\%$	Pass
Infectious Disease NAT HIV 1/2, HEP B/C	HIV 1/2, HEP B/C	Negative
Microbiologic Testing	Not Detected	Not Detected

Stability and Storage

- Follow facility safety protocols when handling and storing cryopreserved cells.
- Store cells in liquid nitrogen vapor phase.
- The product is stable at $< -135^{\circ}\text{C}$ for one year from receipt.
- Short term storage at -80°C may alter the viability of the cells.
- Use samples immediately upon thawing.

Donor Infectious Disease Testing

- HIV 1/2 and Hep B/C testing was performed on maternal blood and/or on a sample of the donated cord blood.
- Viral testing does not guarantee that the donor is virus free. Treat the sample as potentially infectious.

Donor Informed Consent

- Cord blood collection was with consent > approved by an Institutional Review Board and/or in accordance with local, state and national regulatory requirements.

Precautions

- Revvity cannot guarantee the performance of the cells in a researcher's individual assay or culture systems.
- Revvity guarantees that the cells meet the lot specific specifications when assessed immediately after thawing (pre-wash) using flow cytometry methods.

Directions For Use

- To confirm the number, viability and purity of the product, testing must be done immediately post thaw and pre-wash.
- It is recommended that work be performed in a biosafety hood.
 - Wipe the outside of the vial containing the cells with 70% ethanol.
 - Twist the cap slowly to relieve any internal pressure and re-tighten.
 - Thaw cells at 37°C. Do not vortex the cells.
 - Wipe the outside of the vial with 70% ethanol or isopropanol.
 - Determine the volume of the cell suspension.
 - Determine the post thaw cell number, purity and viability based on laboratory standard operating procedures.
 - Wash cells based on laboratories standard operating procedures to prepare cells for downstream applications.
 - Cell loss occurs during wash steps. Do not disturb the cell pellet.

revvity

Revvity, Inc.
77 4th Avenue
Waltham, MA
02451 USA

(800) 762-4000
www.revvity.com

Copyright ©2024, Revvity, Inc. All rights reserved.

FOR RESEARCH ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.