

Research with confidence. Research with Revvity.

Your research deserves a reliable source of cells to support it. Our highly pure and viable stem cells are an ideal choice for a wide range of research and can become an indispensable tool in the lab.

We provide:

- CD34+ cells isolated from human umbilical cord blood
- Exceptional viability (>95%) and high purity (>90%)
- Various cell quantities to help you meet reach your research goals

Working together to support the development of novel applications leading to potential clinical breakthroughs.

CD34+ hematopoietic stem cells from human umbilical cord blood (frozen)



Research ready

Off-the-shelf and ready to use. Let us be your trusted supplier of cells.



Experience matters

Highly qualified and experienced technicians in cryopreservation of cells.



Certificate of analysis

Provided with every sample, giving you the details you need to research with confidence.



Customer support

You and your research outcomes are our top priority. Our responsive team is here to help.

CD34+ hematopoietic stem cells from human umbilical cord blood (frozen)

As a trusted and reliable source of cells, you can plan your experiments with confidence and focus on what really matters – your research outcomes.

Description

- CD34, a transmembrane phosphoglycoprotein, is an established marker of hematopoietic stem and progenitor cells.
- Primary human CD34+ cells are isolated from cord blood using positive immunomagnetic techniques.
- Umbilical cord blood is collected using IRB protocols and/or in accordance with local, state, and federal requirements.
- Cord blood is collected in anticoagulant citrate phosphate dextrose.

Key features

- High purity: > 90% purity by flow cytometry
- Exceptional viability: > 95% cell viability post thaw
- Microbial and infectious disease testing (HIV-1, HIV-2, Hepatitis B, Hepatitis C)
- CD34+ cells frozen in 10% DMSO cryopreservation medium
- Store in liquid nitrogen vapor phase
- Cells stable at < -135°C for one year from receipt
- Certificate of Analysis included

Benefits of working with CD34+ cells

CD34+ cells derived from umbilical cord blood (UCB) offer several advantages compared to those obtained from other sources of CD34+ cells including:

- Non-invasive collection
- More naïve/primitive than other sources
- Greater proliferation

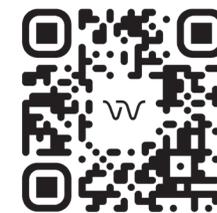
Available cell quantities

- 1 million cells – single donor
Part Number: CBCD34+750-S
- 1 million cells – single donor
Part Number: CBCD34+1M-S
- 2 million cells – single donor
Part Number: CBCD34+2M-S
- 2 million cells – pooled donor
Part Number: CBCD34+2M-P



Potential research applications:

- Molecular genetics
- Angiogenesis
- Regenerative medicine
- Immune diseases
- Gene editing/CRISPR
- Validating expansion technology



revvity