

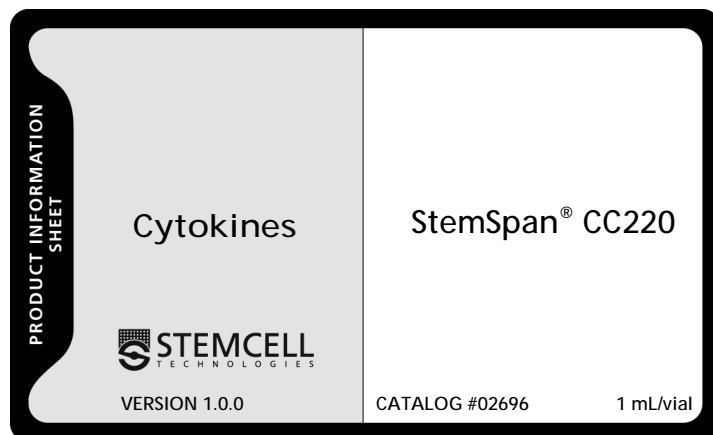
PRODUCT DESCRIPTION

The StemSpan® CC220 cytokine cocktail contains a combination of human recombinant (rh) cytokines formulated to support the expansion of megakaryocyte progenitors and the production of megakaryocytes and platelets in cultures of human hematopoietic cells.

We recommend StemSpan® CC220 in combination with StemSpan® SFEM (Catalog #09600/09650) or StemSpan® H3000 (Catalog #09800/09850) serum-free medium.

StemSpan® CC220 is aseptically manufactured using tightly controlled processes.

Each batch of StemSpan® CC220 is sterility tested. A Certificate of Analysis is available upon request.



FORMULATION

StemSpan® CC220 is provided as a 100X concentrate and contains:

- Recombinant Human Stem Cell Factor (SCF)
- Recombinant Human Thrombopoietin (TPO)
- Recombinant Human Interleukin-6 (IL-6)
- Recombinant Human Interleukin-9 (IL-9)

DIRECTIONS FOR USE

Thaw StemSpan® CC220 at room temperature or at 37°C until just thawed and mix well. If necessary, centrifuge for 30 seconds to remove liquid from cap.

Aseptically transfer StemSpan® CC220 to culture medium at a 100X dilution.

StemSpan® CC220 can be aseptically aliquoted into smaller volumes and stored at -20°C until use.

STABILITY AND STORAGE

StemSpan® CC220 is stable at -20°C until expiry date indicated on label.

StemSpan® CC220 is stable for one month if stored at 4°C under sterile conditions.

Avoid repeated freezing and thawing.