

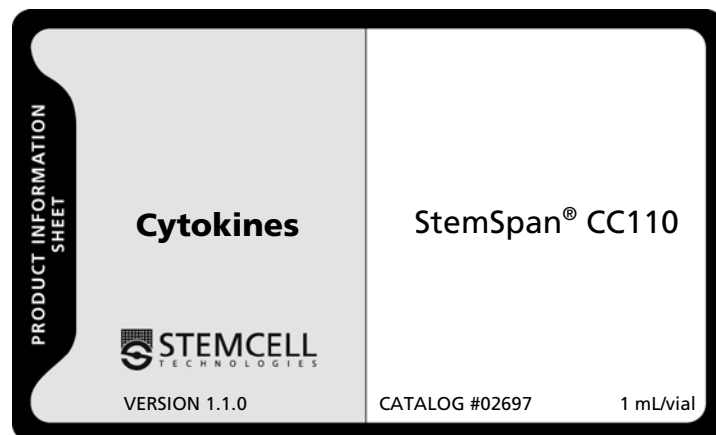
PRODUCT DESCRIPTION

The StemSpan® CC110 cytokine cocktail contains a combination of recombinant human (rh) cytokines formulated to support the proliferation of human hematopoietic progenitors.

We recommend that StemSpan® CC110 be used in combination with StemSpan® SFEM (Catalog #09600/09650) or StemSpan® H3000 (Catalog #09800/09850) serum-free medium.

StemSpan® CC110 is aseptically manufactured using tightly controlled processes.

Each batch of StemSpan® CC110 is sterility tested. A Certificate of Analysis is available upon request.



FORMULATION

StemSpan® CC110 is provided as a 100X concentrate and contains:

- Recombinant Human Flt-3 Ligand (Flt-3-L)
- Recombinant Human Stem Cell Factor (SCF)
- Recombinant Thrombopoietin (TPO)

DIRECTIONS FOR USE

Thaw StemSpan® CC110 at room temperature (15 - 25°C) or at 37°C until just thawed and mix well. If necessary, centrifuge for 30 seconds to remove liquid from cap.

Aseptically transfer StemSpan® CC110 to culture medium at a 100X dilution.

StemSpan® CC110 can be aseptically aliquoted into smaller volumes and stored at -20°C (-25°C to -15°C) until use.

STABILITY AND STORAGE

StemSpan® CC110 is stable at -20°C (-25°C to -15°C) until expiry date indicated on label.

StemSpan® CC110 is stable for one month if stored at 4°C (2 - 8°C) under sterile conditions.

Avoid repeated freezing and thawing.