

## PLTGold<sup>®</sup>

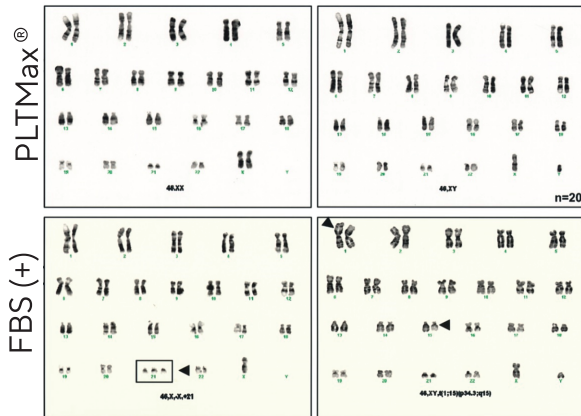
### Human Platelet Lysate

#### Excellent Alternative to FBS

- 사람 혈소판 성분의 FBS 대체제
- Xeno-Free, Heparin-Free
- Donor pool 추출로 Lot-to-lot consistency 확보
- 소량 첨가 (2~5%) 에도 탁월한 세포 증식 효과
- **FDA Type II Master File**
- 40 건 이상의 글로벌 임상시료 생산 레퍼런스 보유
- 품목 구분
  - Research
  - GMP
  - GMP-PI (Pathogen Inactivated)
  - GMP-GI (Gamma Irradiated)



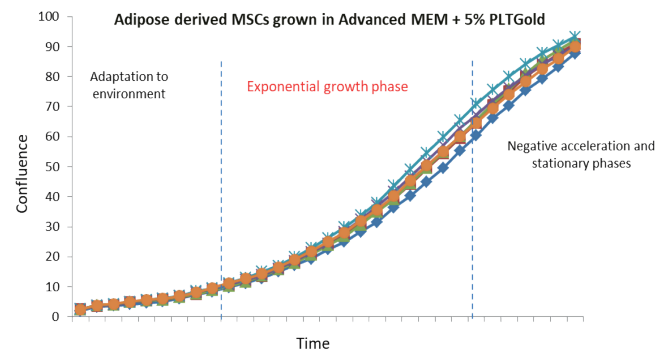
## Early work showed Mill Creek's hPL improved genetic stability over FBS



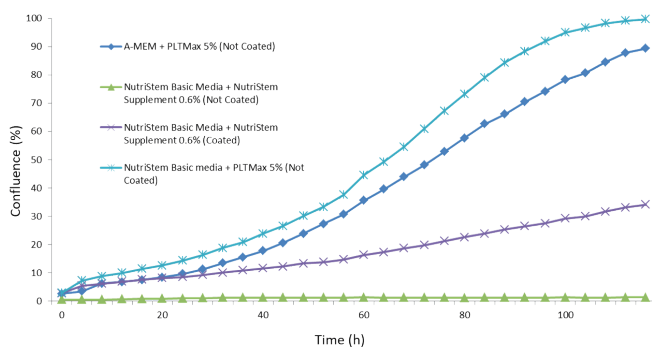
20 adipose derived human MSC were split and cultured for 12 passages in either PLTMax media or FBS. Two FBS cultured samples showed abnormal karyotype. None of the PLTMax samples showed any alterations.

Crespo-Diaz R, et al. Cell Transplant. 2011; 20(6):797-811.

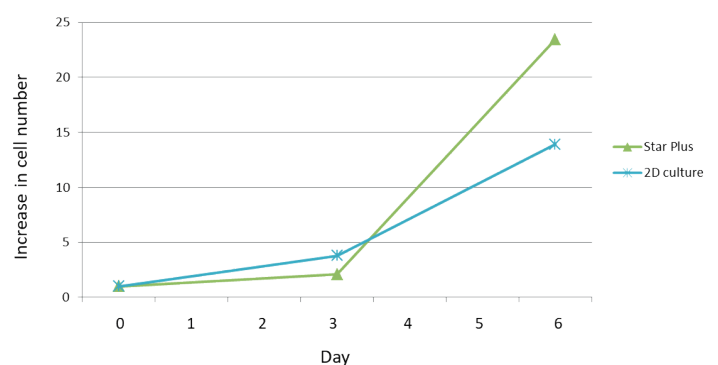
## Lot to lot consistency of PLTGold®



## Early Work Showed PLTMax® enhanced MSC Growth in NutriStem® and eliminated the need for cell attachment coatings



## MSC NutriStem® XF Basal Medium supplemented with PLTGold® and Sartorius microcarriers : Suspension culture



## Regulatory/Quality

- MF on file with the FDA for all Mill Creek clinical grade products
- All products are manufactured under cGMP conditions ISO 9001 (2015) Certification
- Mill Creek has passed regular customer quality audits
- All platelet donors are tested for the required panel of infectious diseases at the time they donate
- Sterility test per USP <71>, Endotoxin per USP <85> and a wide range of other testing is on Product COA
- A panel of growth factors are tested for reference on each lot

## PLTGold® Tips & Tricks

- When the product is received, thaw at 37 °C or 4 °C and aliquot into the sizes you will be using. Freeze what you don't use immediately for later. Too many repeated temperature changes can affect the integrity of the components.
- While we DO NOT recommend filtering PLTGold® or the complete media containing PLTGold®, complete media can be filtered, if necessary.
- We recommend starting with 5% vol/vol and titrating up/down from there to find the sweet spot for each individual application.
- Cell seeding should be performed following the general guidelines for the specific cell type. For Mesenchymal Stem Cells (MSCs), cells are typically plated at approximately  $2 \times 10^3$  –  $5 \times 10^3$  cells per  $\text{cm}^2$ .
- PLTGold® does not contain heparin or require it. DO NOT add heparin.

싸토리우스코리아바이오텍(주)

경기도 성남시 분당구 판교역로 220 쏘리드스페이스 8층

☎ 031. 622. 5700 ✉ marketing@sartorius.co.kr 🌐 www.sartorius.co.kr