

FBS보다 합리적인 선택,
hPL 한 병으로 실험 효율은 UP! 비용은 DOWN!

PLTGold 1병 구매 시, DMEM/RPMI 20병 추가 증정!



Free Gifts!!!

DMEM/ RPMI 20병



x 20



▲ QR 스캔하여 신청하기

PLTGold : Human Platelet Lysate



◀ Datasheet

- 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-GI (Gamma Irradiated)

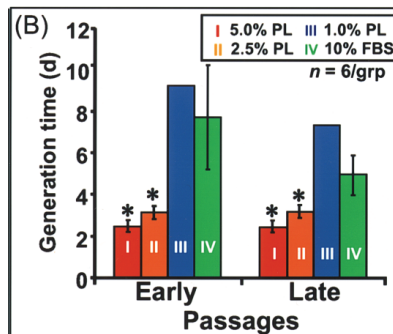
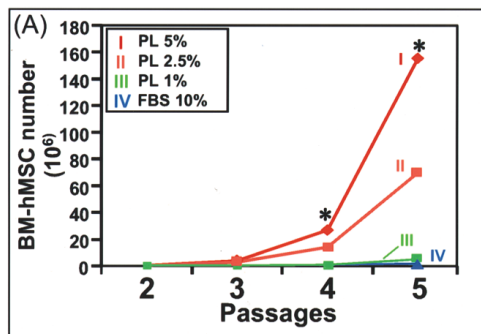
적용 제품

Order No.	Description
PLTGOLD500R	PLTGold Research Grade 500mL
01-050-1A	DMEM, LG, w/o L-Glutamine
01-052-1A	DMEM, HG, with L-Glutamine
01-055-1A	DMEM, HG, w/o L-Glutamine
01-100-1A	RPMI1640 with L-Glutamine
01-104-1A	RPMI1640 w/o L-Glutamine

프로모션 적용 가능 기간 : ~ 2025.12.31

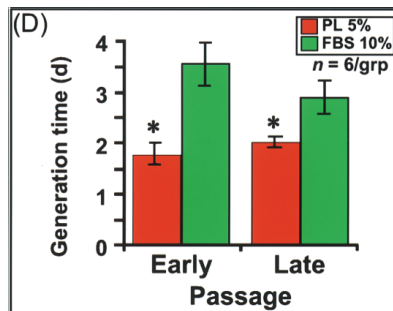
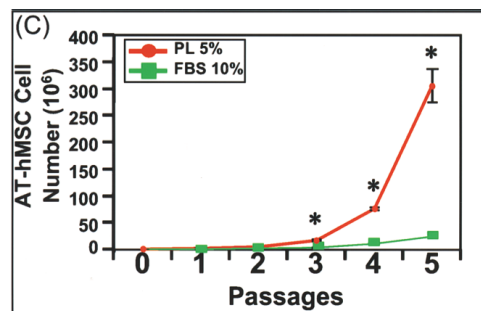
PLTGold® Specifications

Comparison between hPL and FBS of hMSC culture



(A) Bone marrow-derived hMSC (BM-hMSC) cultured in fetal bovine serum (FBS) and different concentrations (1%, 2.5%, and 5%) of platelet lysate (PL) with cumulative quantitation for 5 passages.

(B) Time needed for two population doublings (Generation time) was calculated during early (1-3) and late (3-5) passage for BM-hMSC in the different PL and FBS culture conditions.

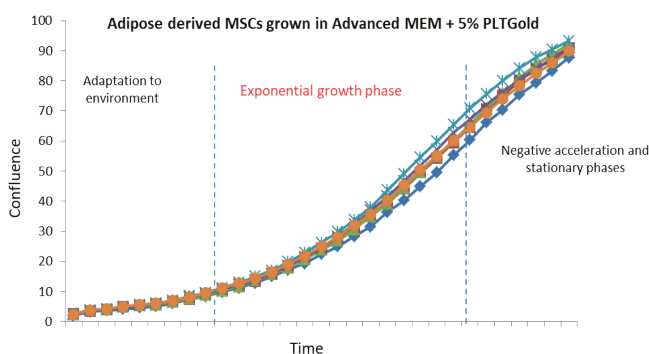


(C) Culture of AT-hMSCs in FBS and 5% GMP-hPL for 5 passages with cumulative quantitation of cellular yield.

(D) Generation time of AT-hMSCs quantified in the GMP-hPL and FBS condition at early (1-3) and late (3-5) passages. *p < 0.01 (A-C) and *p < 0.05 (D).

출처: Crespo-Diaz, Ruben, et al. "Platelet lysate consisting of a natural repair proteome supports human mesenchymal stem cell proliferation and chromosomal stability." Cell transplantation 20.6 (2011): 797-812.

Lot to lot consistency of PLTGold®



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

