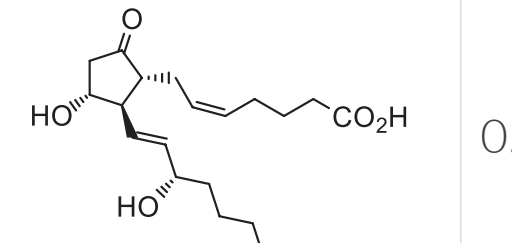
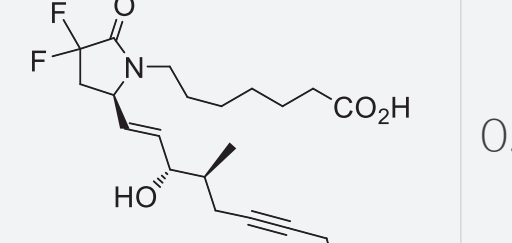


Introduction

In many animals, including humans, PGE₂ causes bone formation with its anabolic actions primarily due to its binding to the GPCR PTGER4 (EP₄).¹ However, systemic PGE₂ administration causes undesired side effects such as pain, lethargy, and diarrhea due to its binding to the EP₁ and EP₃ receptors.² The current bone anabolic therapeutic agents are primarily protein based, requiring appropriate storage conditions and having relatively short shelf lives. They are also expensive, and the morphogenic activity associated with many of them raises safety concerns and limits their potential use in orthopedic and other local applications. To alleviate these issues in a potential therapeutic, we developed a series of novel difluorolactam EP₄ receptor agonists.³ The work presented here is from the ongoing characterization of our lead compound, KMN-159, which shows high selectivity for EP₄ binding, excellent potency for EP₄ activation, potent stimulation of osteoblastic differentiation in unfractionated bone marrow cells (BMCs), and remarkable chemical stability.

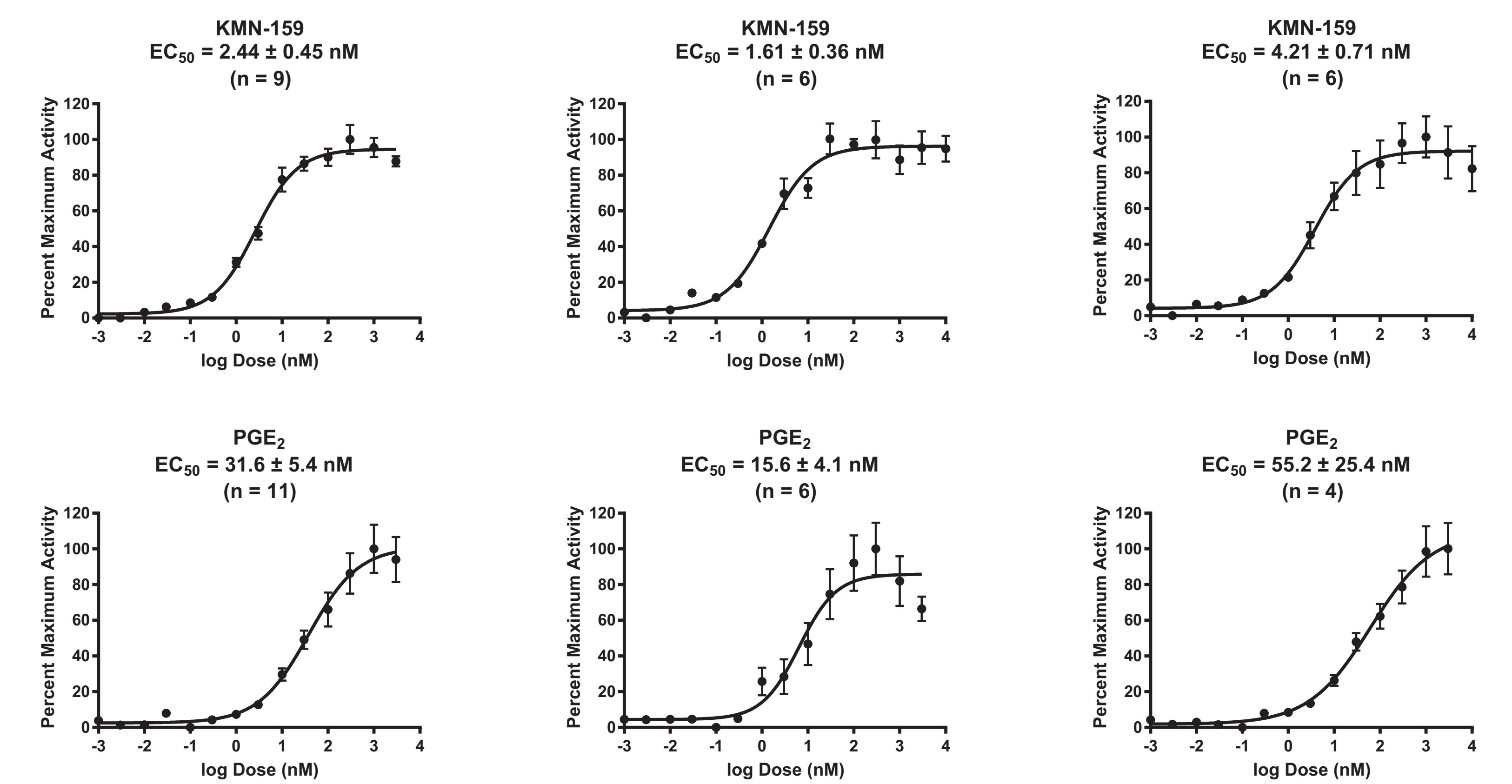
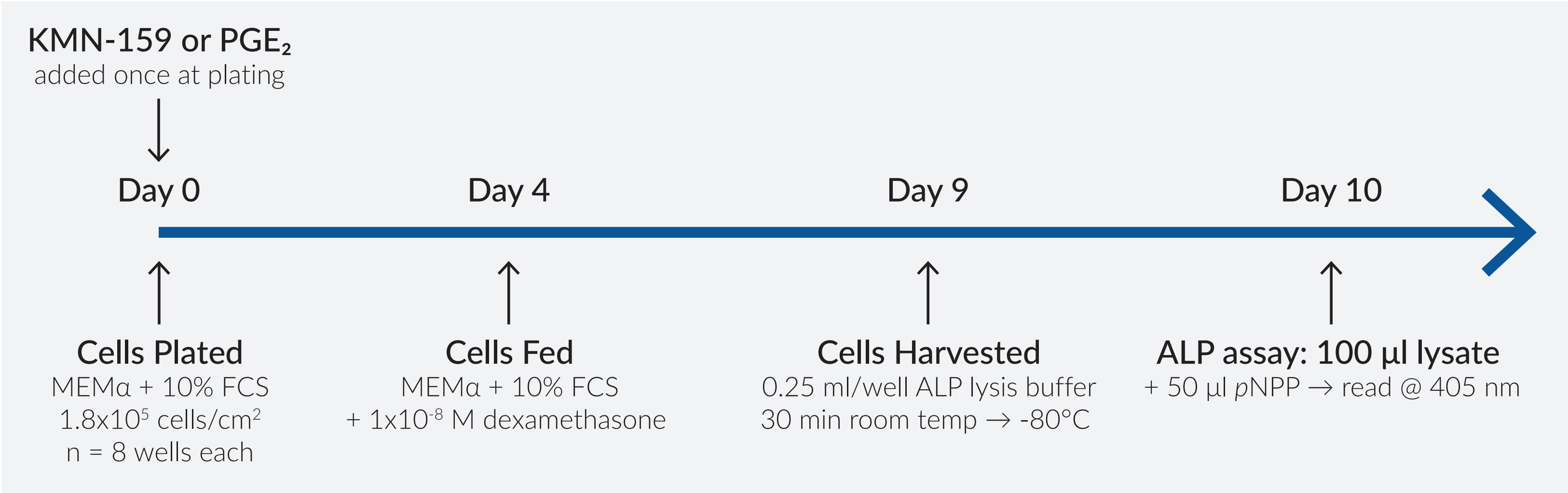
KMN-159 is a Highly Selective and Potent Activator of the Rat EP₄ Receptor

While the natural ligand PGE₂ is a potent activator of both the rat EP₂ and EP₄ receptors, KMN-159, the lead compound in our novel difluorolactam series, demonstrates potency for binding, stimulation of cAMP production, and activation of a CRE-SEAP reporter that is equivalent to PGE₂ but does so in a clearly EP₄-selective manner.

Compound	Structure	rat EP ₄ receptor binding K _i (nM)	rat EP ₄ receptor cAMP formation EC ₅₀ (nM)	rat EP ₄ receptor CRE activation EC ₅₀ (nM)	rat EP ₄ receptor cAMP formation EC ₅₀ (nM)	rat EP ₄ receptor CRE activation EC ₅₀ (nM)
PGE ₂		0.19	88.4 ± 4.2 (n = 5)	2.8 ± 0.7 (n = 9)	0.48 ± 0.09 (n = 5)	0.03 ± 0.002 (n = 56)
KMN-159		0.24	>10,000 (n = 3)	>10,000 (n = 3)	0.50 ± 0.08 (n = 4)	0.03 ± 0.003 (n = 7)

Binding cAMP measurement and CRE activation assays were all performed as in Barrett *et al.*, *J. Med. Chem.*, 2019.³

KMN-159 Stimulates Osteoblastic Differentiation in Young and Old Female and Young Male Rat BMCs



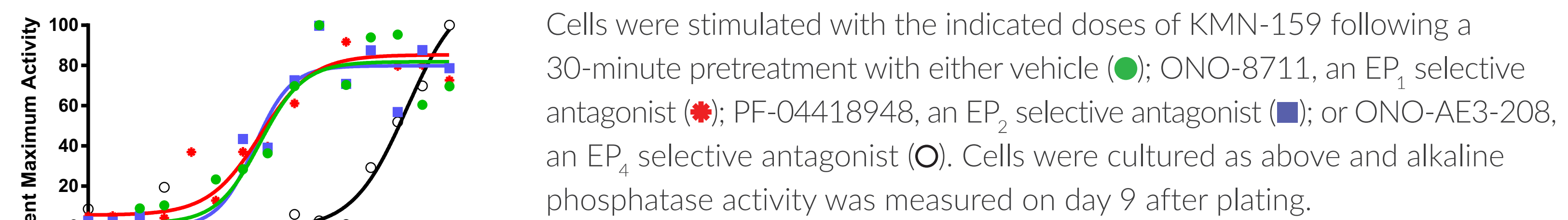
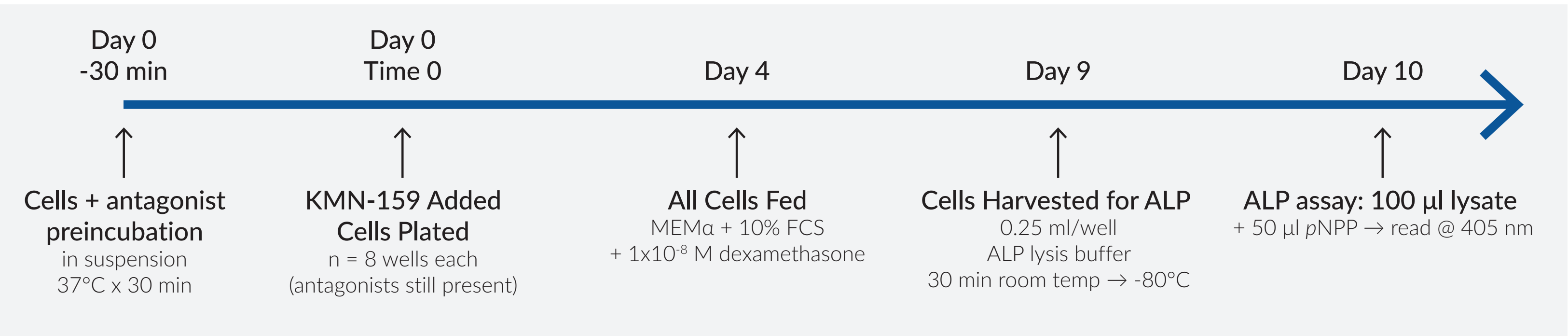
Young Female (<6 months)

Old Female (>12 months)

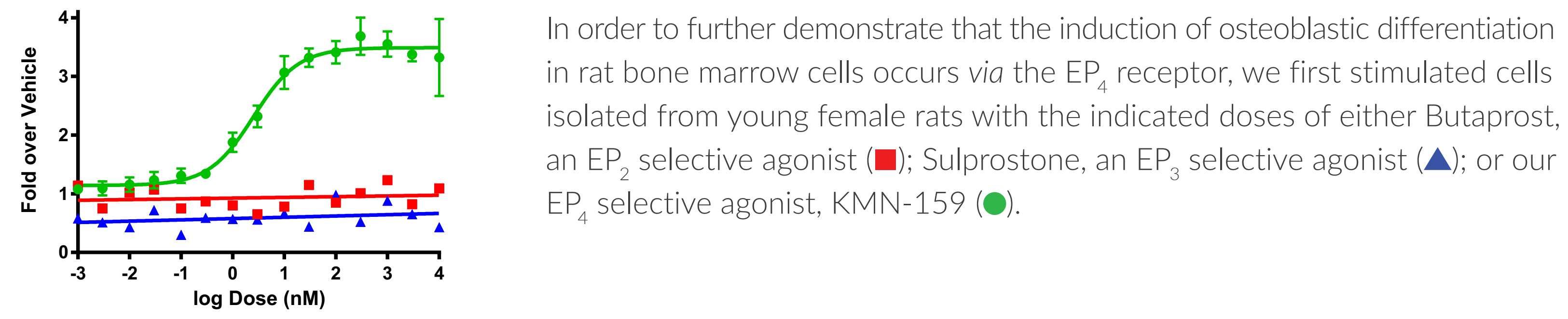
Young Male (<6 months)

Stimulation of Osteoblastic Differentiation in Rat Bone Marrow Cells Occurs via EP₄ Activation

These data demonstrate that KMN-159 stimulates osteogenesis *via* activation of the EP₄ receptor; specific agonists for the related EP₂ and EP₃ receptors do not stimulate osteogenesis and only an EP₄-selective antagonist blocks the osteogenesis induced by KMN-159.

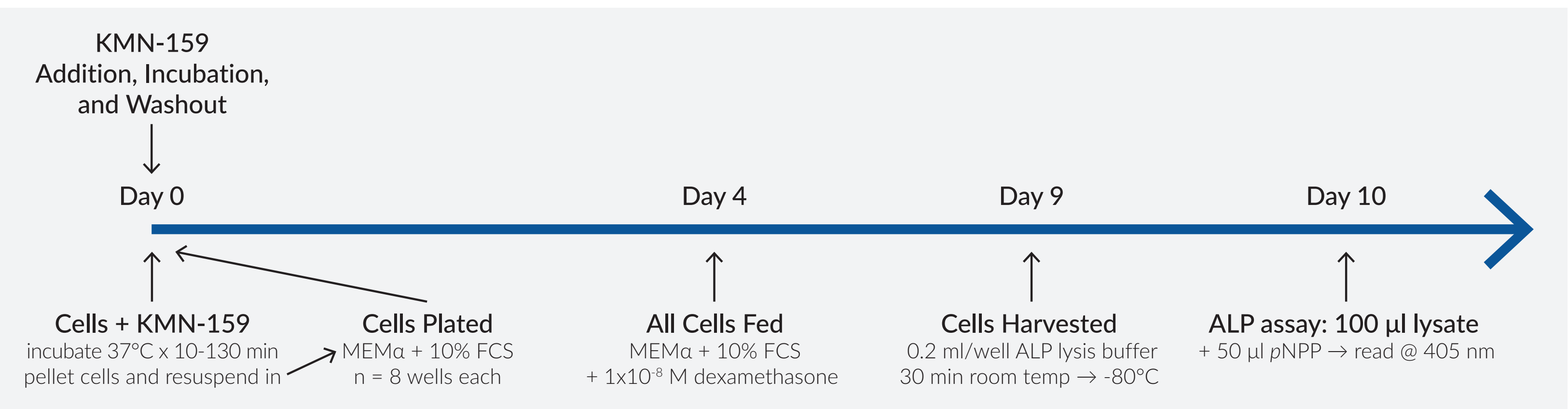


Cells were stimulated with the indicated doses of KMN-159 following a 30-minute pretreatment with either vehicle (●); ONO-8711, an EP₄ selective antagonist (●); PF-04418948, an EP₂ selective antagonist (■); or ONO-AE3-208, an EP₃ selective antagonist (○). Cells were cultured as above and alkaline phosphatase activity was measured on day 9 after plating.



In order to further demonstrate that the induction of osteoblastic differentiation in rat bone marrow cells occurs *via* the EP₄ receptor, we first stimulated cells isolated from young female rats with the indicated doses of either Butaprost, an EP₂ selective agonist (■); Sulprostone, an EP₃ selective agonist (▲); or our EP₄ selective agonist, KMN-159 (●).

Commitment to Osteogenic Differentiation Requires Only a Short Exposure to KMN-159

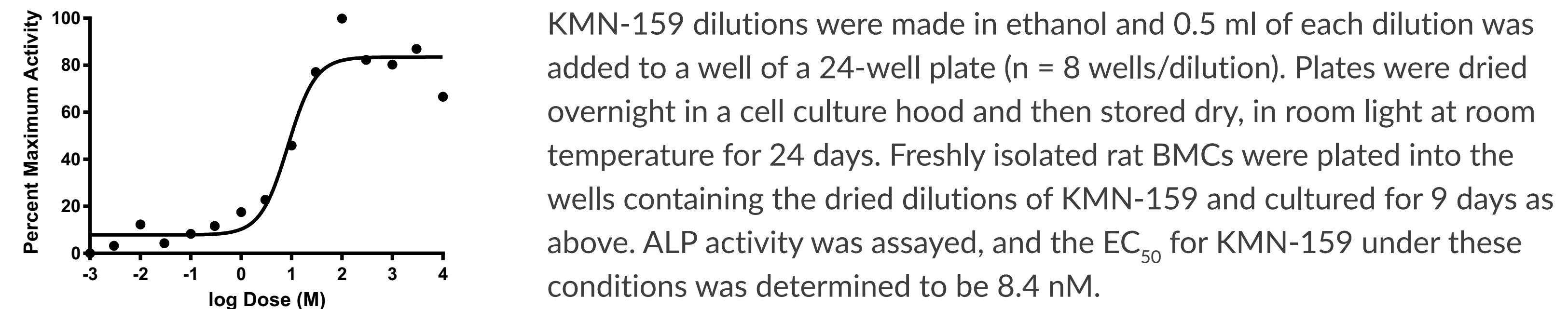


Alkaline Phosphatase (Day 9)

Mineralization (Day 21)

These data show that rat BMCs exposed to KMN-159 for as little as 10 minutes are committed to induction of alkaline phosphatase measured 9 days later. While only 30- and 60-minute exposures of the cells to KMN-159 were tested for mineralization, it is clear that these short exposure times lead to the commitment of stem cells within the bone marrow to the complete program of osteogenesis. This is demonstrated by increased mineralization with 1,000 nM KMN-159 treatment (right half of each plate) as compared to vehicle treatment (left half of each plate). This is consistent with the typical transient activation of a GPCR such as EP₄.

KMN-159 Stimulates ALP Activity Even After Dry Plate Storage at Room Temperature

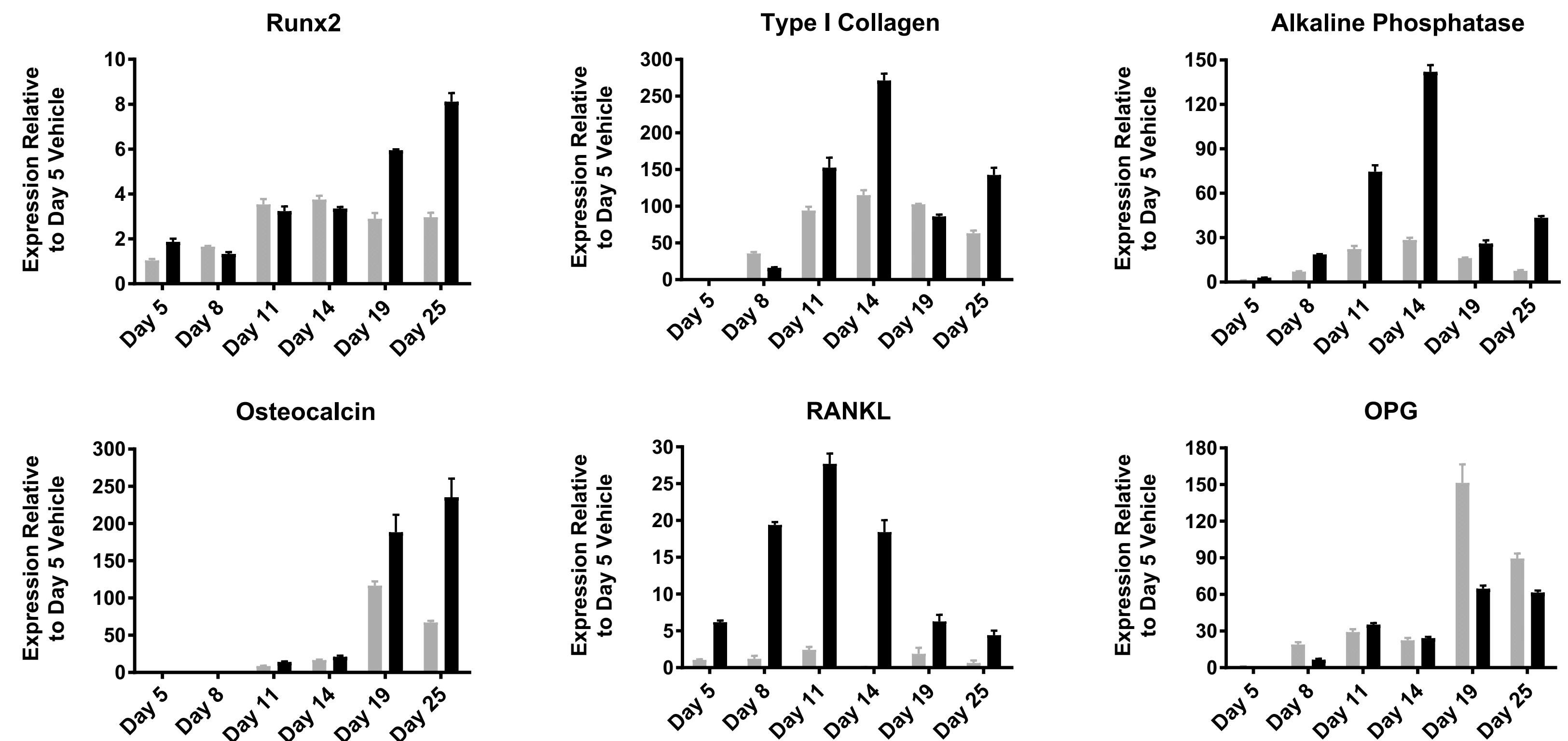


KMN-159 dilutions were made in ethanol and 0.5 ml of each dilution was added to a well of a 24-well plate (n = 8 wells/dilution). Plates were dried overnight in a cell culture hood and then stored dry, in room light at room temperature for 24 days. Freshly isolated rat BMCs were plated into the wells containing the dried dilutions of KMN-159 and cultured for 9 days as above. ALP activity was assayed, and the EC₅₀ for KMN-159 under these conditions was determined to be 8.4 nM.

KMN-159 was active in stimulating ALP enzyme activity in BMCs with an EC₅₀ value equivalent to that seen in cells treated directly with compound. This demonstrates that the compound is stable at room temperature and retains its potency for at least 24 days.

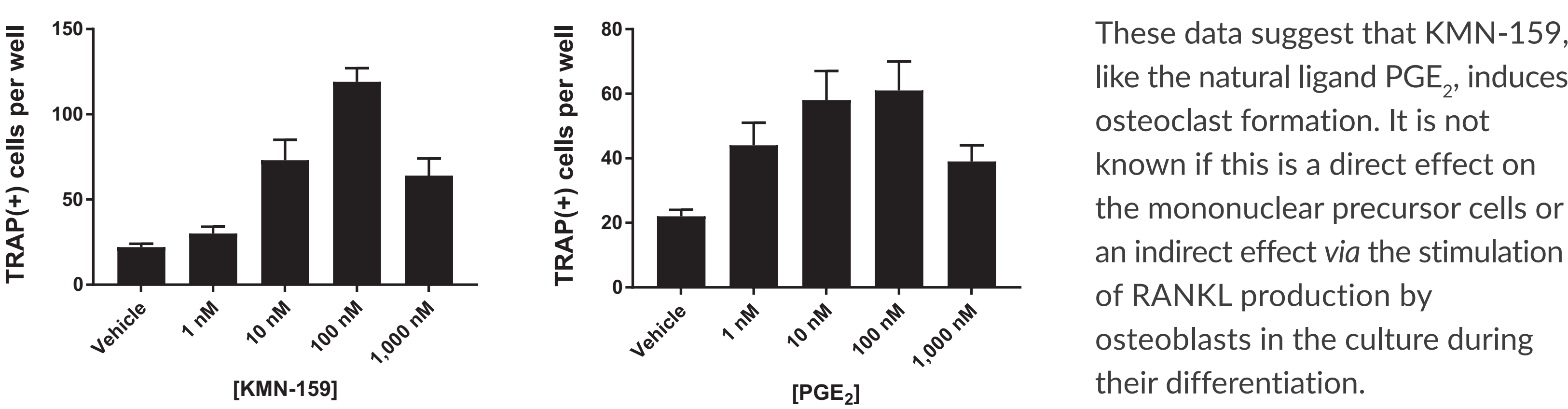
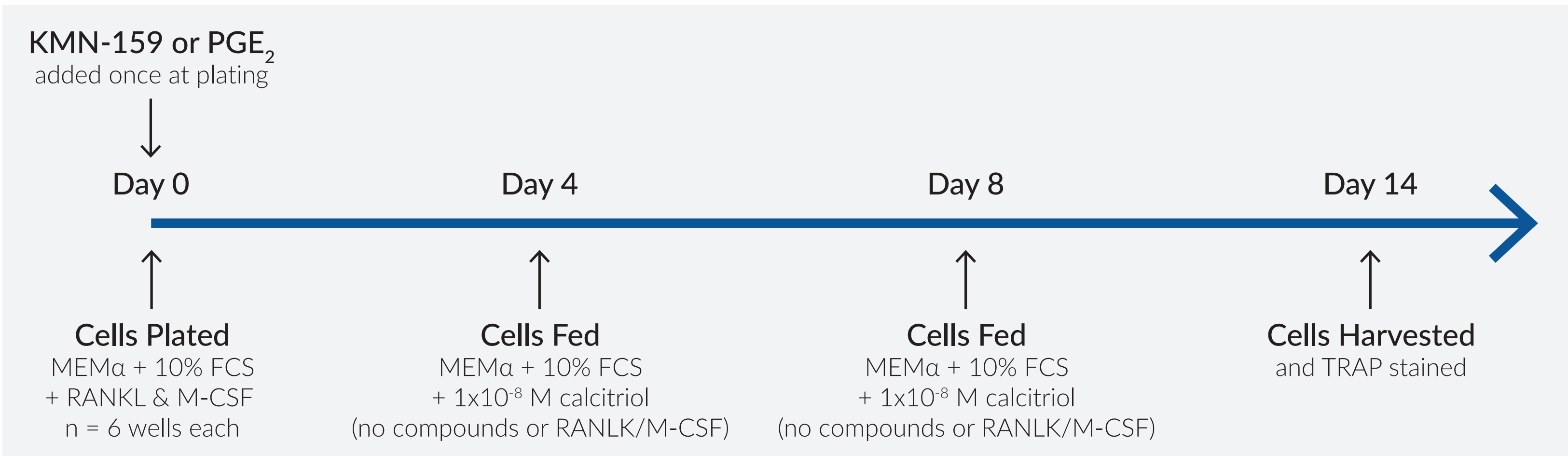
KMN-159 Stimulates Expression of Osteoblast Phenotype Marker Genes

Phenotype marker genes were expressed in a manner consistent with the stimulation of osteoblastic differentiation by KMN-159. RANKL expression was also stimulated, followed by a later increase in OPG expression, consistent with an observed increase in TRAP-positive cells by KMN-159 (see below).

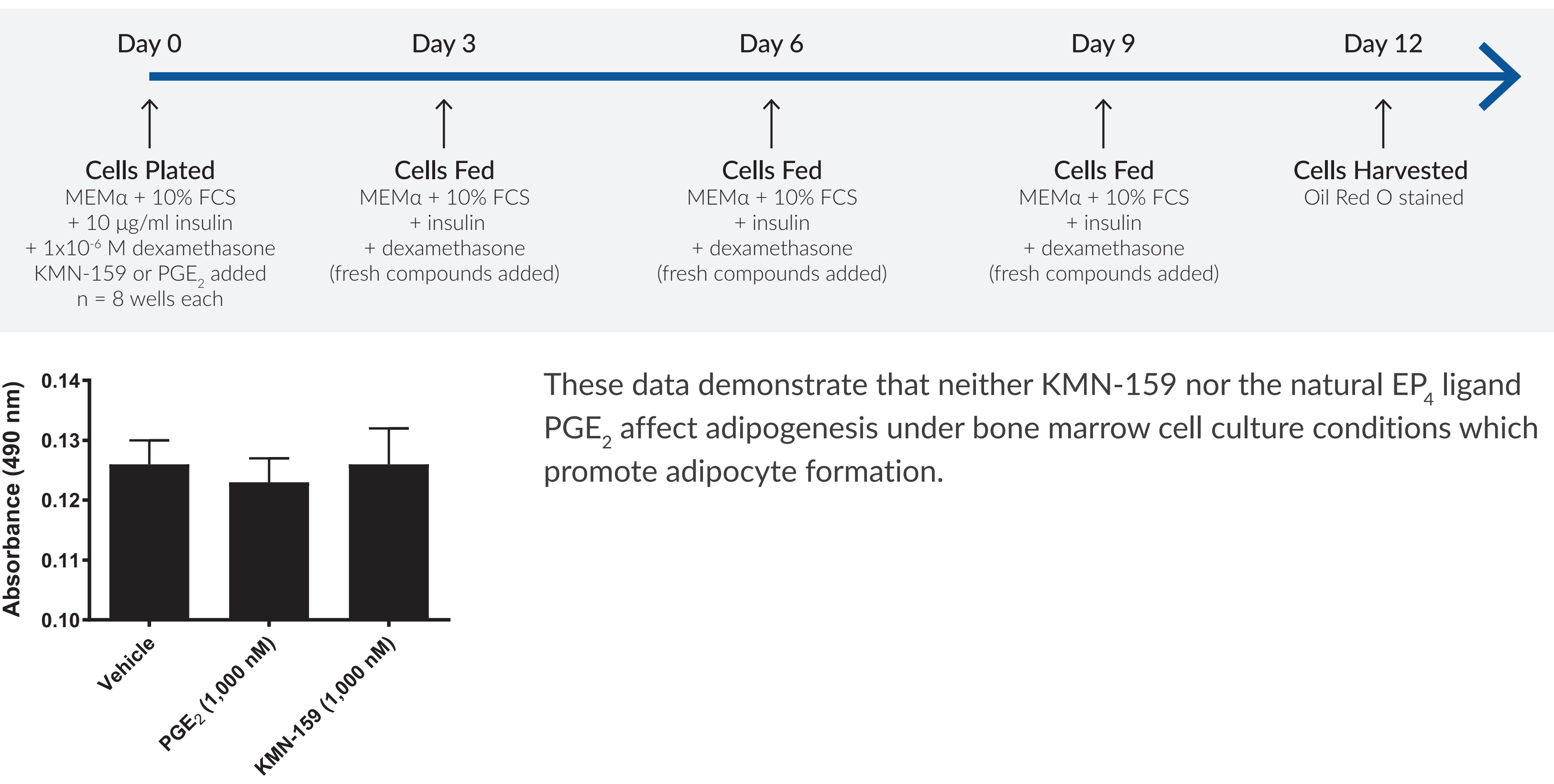


Rat BMCs were treated with 1,000 nM KMN-159 or ethanol vehicle and plated in 12-well dishes as above. Cells were fed every 3 days with osteogenic medium beginning on day 7. RNA was prepared on the indicated days and used as a template for RT-qPCR using GAPDH as a reference gene. (vehicle = ■ KMN-159 = ■)

KMN-159 Stimulates Osteoclastogenesis but not Adipogenesis



These data suggest that KMN-159, like the natural ligand PGE₂, induces osteoclast formation. It is not known if this is a direct effect on the mononuclear precursor cells or an indirect effect *via* the stimulation of RANKL production by osteoblasts in the culture during their differentiation.



These data demonstrate that neither KMN-159 nor the natural EP₄ ligand PGE₂ affect adipogenesis under bone marrow cell culture conditions which promote adipocyte formation.

Conclusions

- KMN-159, a novel difluorolactam EP₄ receptor agonist, specifically binds to and activates the rat EP₄ receptor with potency similar to the natural ligand PGE₂.
- KMN-159 stimulates osteoblastic differentiation in BMCs regardless of the age or sex of the donor rats.
- When dried on the surface of cell culture wells and stored at room temperature for at least 24 days, KMN-159 retains the potency and efficacy equivalent to that seen when cells are treated with freshly prepared stocks.
- Consistent with the EP₄ receptor being a GPCR, BMCs need to be exposed to KMN-159 for less than 30 minutes to fully initiate osteogenesis as measured by ALP and mineralization.
- The osteoblastic differentiation induced by KMN-159 occurs solely through stimulation of the EP₄ receptor as demonstrated by receptor-specific agonist and antagonist studies.
- Osteoblast phenotype marker genes including Runx2, type I collagen, ALP, and OC are expressed at higher levels in the KMN-159 treated cells than in the vehicle treated controls.
- The expression of RANKL is increased by KMN-159.
- KMN-159 and PGE₂ both increased the number of osteoclasts but not adipocytes.

Significance and Future Directions

- The work presented here used whole rat bone marrow to explore the action of KMN-159. This complex mixture of cells is similar to the milieu in which the compound would act as a therapeutic agent for fracture therapy.
- Complete stimulation of osteogenic differentiation in bone marrow cells by a single dose of KMN-159 with an exposure time of less than 30 minutes suggests that a combination of this compound with a surgically appropriate matrix will allow a single clinical application to be performed. This will provide advantages over current therapies including long-term shelf stability and a greatly reduced cost of goods.
- The combination of potency of KMN-159 and its selectivity for EP₄, as shown by several measures, should decrease the risk of off-target activity following release from a matrix in a local administration modality.
- Application of KMN-159 or similar compounds in a local manner to increase the rate of bone formation should have many clinical uses including the augmentation of bone mass prior to implant, spinal fusion, repair of non-union fractures, bridging long bone critical defects, coating implants, and craniofacial repair.

Acknowledgements

- The authors would like to acknowledge the continuing support by the staff members of Cayman Chemical and Ramapo College of New Jersey.
- This work was supported by a grant from the Ramapo College Foundation, the Cayman Intramural Research Project, and the TAS Research Grants Program at Ramapo College.

References
1. Pagliaro, J., Leonidou, A., Ar-Saltany, M., et al. *Curr. Mol. Pharmacol.* 5(2), 174-181 (2012).
2. Li, M., Thompson, D.D., and Paraskas, V.M. *Int. Orthop.* 33(6), 767-772 (2009).
3. Barrett, S.D., Hsiao, M.C., Korman, J.B., et al. *J. Med. Chem.* 62(9), 4751-4761 (2019).