

Novel Potent Lactam Acetylene EP₄ Agonists Stimulate Alkaline Phosphatase Production and Differentiation in Bone Marrow Cells

Stephen D. Barrett,*¹ Fred L. Ciske,¹ Gregory W. Endres,¹ Thomas A. Owen,² Adam Uzieblo,¹ Bradlee D. Germain,¹ Andrei Kornilov,¹ Joseph M. Colombo,¹ James B. Kramer,¹ Liyue Huang,¹ Jeffrey K. Johnson,¹ James P. O'Malley²

¹ Cayman Chemical Company, Inc. Ann Arbor, MI

² Myometrics, LLC New London, CT



► Abstract

The prostaglandin E G-protein coupled receptors (GPCR) subtypes 2 and 4 (EP₂ and EP₄ receptors, respectively) are stimulated by interaction with two eicosanoid products of the cyclooxygenase-catalyzed pathway, prostaglandin E₁ (PGE₁) and prostaglandin E₂ (PGE₂), and signal by increasing intracellular cAMP levels.¹ Expression of EP₂ and EP₄ receptors in human bone and their roles in differentiation, growth, and remodeling have been previously investigated.² Although the presence and function of the EP₂ receptor remains ambiguous, the clear significance of the EP₄ receptor in bone anabolism has been demonstrated with both systemic and locally-administered selective EP₄ agonists. Most osteoporosis programs focused on generating agonists with high systemic exposure and selectivity for EP₂/EP₄ versus EP₁ and EP₃. Substituted γ -lactam (pyrrolidinone) derivatives that structurally mimic the carbocyclic prostaglandin scaffold were found to possess notable selectivity for EP₄ over the other EP receptor subtypes, especially EP₁. As the osteoporosis development programs were failing for safety reasons associated with high systemic exposures, attention shifted to localized delivery of anabolic agents to bone, resulting in the approval of the costly recombinant human bone morphogenetic protein 2 (BMP-2) for orthopedic indications. We have prepared novel γ -lactam derivatives designed to maximize selectivity for EP₄ versus EP₁, EP₃, and other prostanoid receptors and to minimize systemic exposure. These compounds were first screened in EP₄ and EP₂ transfected cells by measuring their effect on secreted alkaline phosphatase (SEAP) stimulation. Compounds of interest were then tested for their ability to stimulate osteoblastic differentiation of primary rat bone marrow cells.

► IN VITRO SCREENING ASSAYS

RECEPTOR BINDING ASSAYS

EP₁-EP₄: [³H]PGE₂ displacement/competition assay of human prostanoid EP₁₋₄ receptor in transfected HEK-293 cells (Cerep, Catalog Reference Nos. 0440, 1955, 2774, 0441, respectively)

FP: [³H]PGF_{2a} displacement/competition assay of human prostanoid FP receptor in transfected HEK-293 cells (Cerep, Catalog Reference No. 1979)

DP₁: [³H]PGD₂ displacement/competition assay of human prostanoid DP₁ receptor in transfected 1321N1 cells (Cerep, Catalog Reference No. 2517)

IP: [³H]iloprost displacement/competition assay of human prostanoid IP receptor in transfected HEK-293 cells (Cerep, Catalog Reference No. 2230)

FUNCTIONAL CELL ASSAYS (STEP Plate Format)

Both SEAP activity and cAMP assays for EP₂ or EP₄ agonists were performed on EP₂ or EP₄ STEP (Surface Transfection and Expression Protocol) plates (from Cayman) which are coated with rat EP₂ or EP₄ receptor and secreted alkaline phosphatase (SEAP) -cyclic AMP response element (CRE) reporter constructs. Cells grown on the STEP complex will express EP₂ or EP₄ at the cell surface. Binding of agonists to EP₂ or EP₄ initiates a signal transduction cascade results in a transient increase in cAMP and an prolonged increase in expression of SEAP which is secreted into the cell culture media. cAMP levels were then measured by ELISA and SEAP activity was measured with a luminescence-based alkaline phosphatase substrate.

EP₂ and EP₄ Transfected HEK-293 Cell Assay (SEAP)

1. Cells were seeded on an EP₂ or EP₄ STEP plate at a density of 40,000 – 80,000 cells/well in 200 μ L of reduced serum medium containing 0.5% FBS and incubated at 37 °C incubator with 5% CO₂ for 16-18 hour.

2. Culture media was aspirated and 200 μ L of culture medium containing test compounds was added to wells. For each compound, an 8-point dose response curve (DRC) ranging from 10 μ M to 0.001 pM was performed in triplicate. PGE₂ DRCs were run in parallel in all experiments (concentrations from 0-6,000 pM for EP₄ and 0-10,000 nM for EP₂)

3. After 6-8 hours of stimulation, 10 μ L of media was transferred to a corresponding well of a 96-well solid black plate. The plate was heated at 65 °C for 30 minutes to inactivate endogenous alkaline phosphatase.

4. Luminescence-based alkaline phosphatase substrate (Michigan Diagnostics, LLC, Cat#SAP450101) was added to each well and SEAP activity was measured by reading the luminescent signal.

5. The EC₅₀ values for PGE₂ and each test compound were calculated using GraphPad Prism 5.

EP₂ and EP₄ Transfected HEK-293 Cell Assay (cAMP)

- As above.
- As above, except the culture medium contained 500 μ M IBMX (an inhibitor of cAMP phosphodiesterase).
- Cells were incubated at 37 °C for 30 minutes.
- The plate was centrifuged at 1,000 x rpm for 10 minutes and the supernatant aspirated.
- EIA assay buffer (100 μ L) was added to each well and cells were lysed by freezing at -80 °C freezer for 1 hr followed by completely thawing at room temperature.
- The plate was centrifuged at 1,000 x rpm for 10 minutes. The supernatant was assayed for cAMP by ELISA (Cayman Chemical item 581001).
- The data were analyzed using GraphPad Prism 5 as above.

TP TRANSFECTED HEK-293 CELL ASSAY (intracellular [Ca²⁺])

Cerep Catalog Reference No. 2059

► TABLE 1 ω -Chain SAR

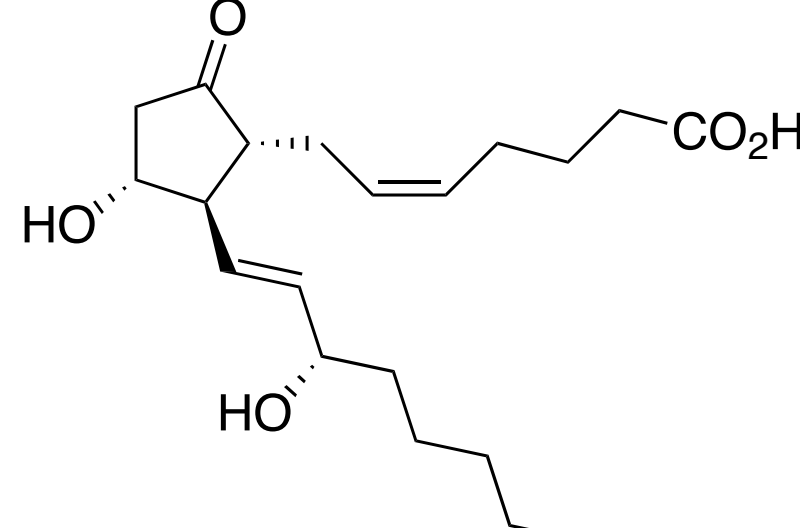
		PGE ₂ Displacement Radioligand Binding EP ₄ IC ₅₀ (nM)	HEK-293 Cells SEAP Stimulation EP ₄ EC ₅₀ (nM)	Bone Marrow Cells EC ₅₀ (nM)
PGE ₁			0.22 ± 0.04 (cAMP)	
KMN-182		2.0	0.71 (cAMP)	
KMN-42			23.5	
KMN-100			4.73	
KMN-64			371	
KMN-32		3.9	0.57	
KMN-80		3.0	0.19 ± 0.07	
KMN-285		54	0.65	
KMN-293		>100	0.33	
KMN-33			4.05	
KMN-79		6.9	1.14	
KMN-287		4.8	3.96	
KMN-295		>100	7.72	

► TABLE 2 α -Chain SAR

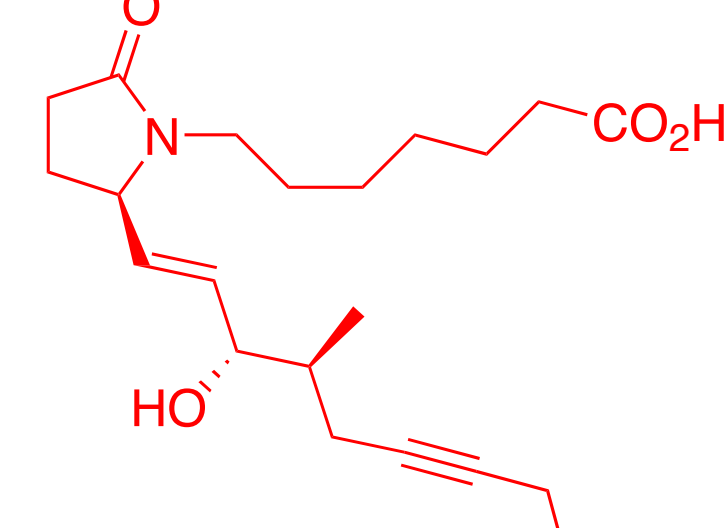
		PGE ₂ Displacement Radioligand Binding EP ₄ IC ₅₀ (nM)	HEK-293 Cells SEAP Stimulation EP ₄ EC ₅₀ (nM)	Bone Marrow Cells EC ₅₀ (nM)
PGE ₂		0.38 ± 0.07	0.05 ± 0.03	
KMN-151			2.57	
KMN-123		0.67	0.31	
KMN-212		6.1	1.54	

► TABLE 3

PGE₂, KMN-80, Compound X and Compound Y prostanoid profiles



PGE₂



KMN-80

Compound X

Compound Y

Radioligand Binding IC₅₀ (nM)

	EP ₁	EP ₂	EP ₃	EP ₄	FP	DP ₁	IP
PGE ₂	1.4	2.6	1	2.0	199 ± 12 ³	307 ± 106 ³	>10,000 ³
KMN-80	>10,000	>10,000	1,400	3.0	>10,000	>10,000	>10,000
Compound X	>10,000	9,000	430	1.3	>10,000	>10,000	>10,000
Compound Y	>10,000	120	400	0.74	>10,000	>10,000	>10,000

Functional EC ₅₀ (nM)			
	EP ₂	EP ₄	TP
PGE ₂	59 ± 17	0.05 ± 0.03	>20,000 (binding K) ³
KMN-80	>1,000	0.19 ± 0.07	>10,000
Compound X	>1,000	0.04 ± 0.04	~10,000
Compound Y	394	0.01	>10,000

Bone Marrow Cells		
EC ₅₀ (nM)		
PGE ₂	27	Vehicle
KMN-000080	20	
Compound X	3	Compound Y (30 nM)
Compound Y	0.29	

► PROGRAM GOALS

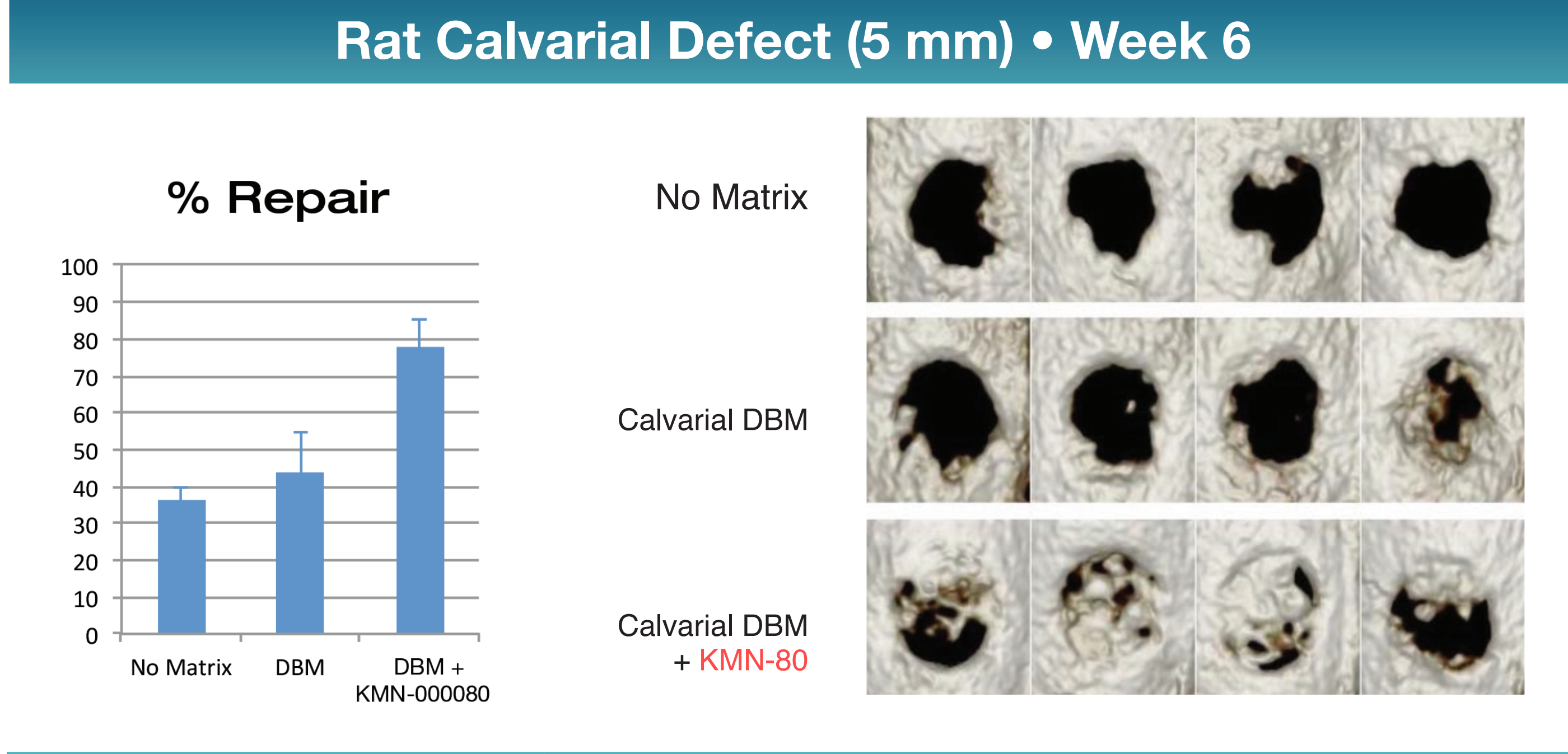
- Novel, small molecule, potent EP₄ receptor agonist
- Selectivity for the EP₄ receptor versus other prostanoid receptors
- Suitable physical property and pharmacokinetic profile for local administration
- Efficacy at tolerated doses in combination with delivery matrix

► PROGRAM ACHIEVEMENTS

- KMN-80, Compound X, and Compound Y are potent in EP₄ receptor binding and functional/cellular assays.
- Selective for the EP₄ receptor against all other prostanoid receptors
- Soluble in local-administration matrices, and Compound Y is highly cleared (low systemic exposure at efficacious doses)
- Efficacious in key animal models; Compound Y is Ames negative

► FIGURE 1

Rat Calvarial Defect treated with KMN-80 in demineralized bone matrix (DBM)



RAT BONE MARROW CELL ASSAY

- Extrude bone marrow from both tibia and femora of intact female rats between 6 and 12 weeks of age (Harlan) into complete cell culture medium (MEMa) supplemented with 15% fetal calf serum and antibiotics), filter through a 100 μ m mesh filter, and count.
- Plate cells into either 24-well (for alkaline phosphatase assay) or 6-well (for nodule formation) dishes at 1.78x10⁵ cells/cm² in complete medium.
- Culture cells (37 °C, 5% CO₂, humidified) for seven days with one media change on day four (50% media withdrawn and replaced with an equal volume of fresh media containing 2x10⁻⁸ M dexamethasone (1x10⁻⁶ M final).
- For the alkaline phosphatase assay, harvest cells at day 7 by incubation in alkaline phosphatase lysis buffer and store at -80 °C. Quantitate alkaline phosphatase activity by incubating 100 μ L cell lysate with 50 μ L pNPP substrate and reading product formation at 405 nm. Determine EC₅₀ values were from the data using GraphPad Prism.
- For nodule formation, feed cells every 3 days beginning at day 7 with MEMa supplemented with 10% FCS, antibiotics, 50 mg/ml ascorbic acid, 2 mM inorganic phosphate, and 1x10⁻⁸ M dexamethasone. At the last feeding on day 18, supplement the medium with calcein (1 mg/ml). Fix the cells in 10% buffered formalin on day 21 and photograph under UV light.

RAT CALVARIA IN VIVO ASSAY

Will be described in more detail in Poster [MO0405] at Poster Session III on Monday, October 7.

REFERENCES

- Fujino, H. and Regan, J., *Trends in Pharmacological Sciences*, **2003**, 24(7), 335-340; Hoshino, T. *et al.*, *J. Biol. Chem.*, **2003**, 278(15), 12752-12758; Takahashi, S. *et al.*, *Biochem. Pharmacol.*, **1999**, 58(12), 1997-2002; Quiroga, J. *et al.*, *Pharmacol. Ther.*, **1993**, 58(1), 67-91.
- Iwaniec, U. *et al.*, *Osteoporosis International*, **2007**, 18(3), 351-362; Aguirre, J. *et al.*, *J. Bone and Min. Res.*, **2007**, 22(6), 877-888; Yoshida, K. *et al.*, *Proc. Natl. Acad. Sci. USA*, **2002**, 99 (7), 4580-4585. Hayashi, K. *et al.*, *J. Bone Joint Surg. Br.*, **2005**, 87-B (8), 1150-6.
- Young, R. *et al.*, *Heterocycles*, **2004**, 64, 437-445. Receptor binding assays were performed using membranes from HEK-293ebna cells recombinantly expressing the corresponding human prostanoid cDNAs.