

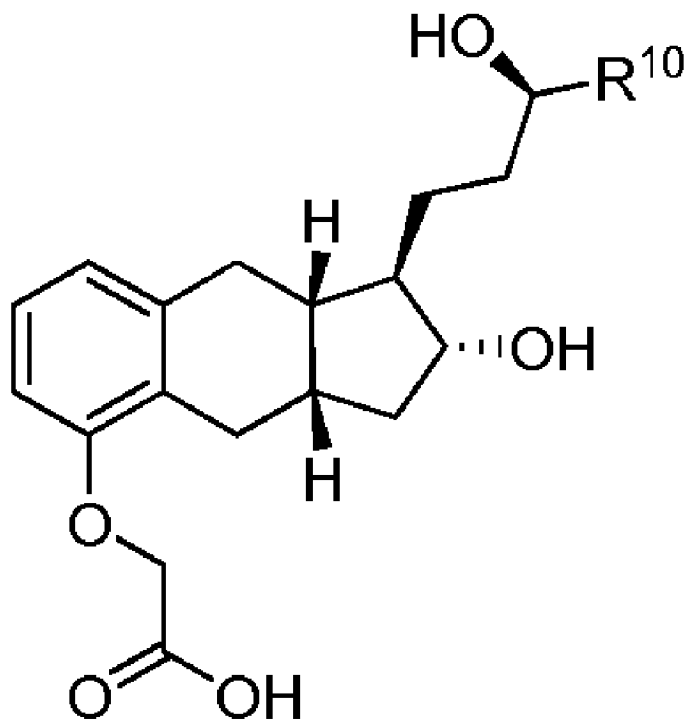


(86) **Date de dépôt PCT/PCT Filing Date:** 2013/12/06
(87) **Date publication PCT/PCT Publication Date:** 2014/06/12
(85) **Entrée phase nationale/National Entry:** 2015/06/01
(86) **N° demande PCT/PCT Application No.:** US 2013/073474
(87) **N° publication PCT/PCT Publication No.:** 2014/089385
(30) **Priorités/Priorities:** 2012/12/07 (US61/734,672);
2013/03/12 (US61/777,882)

(51) **Cl.Int./Int.Cl. C07C 205/57** (2006.01),
C07C 51/347 (2006.01), **C07D 301/32** (2006.01),
C07F 7/18 (2006.01)
(71) **Demandeur/Applicant:**
CAYMAN CHEMICAL COMPANY INCORPORATED, US
(72) **Inventeurs/Inventors:**
HERING, KIRK WILLIAM, US;
CHAMBOURNIER, GILLES, US;
ENDRES, GREGORY WILLIAM, US;
FEDIJ, VICTOR, US;
KRELL, THOMAS JAMES, II, US;
MAHMOUD, HUSSEIN MAHMOUD, US
(74) **Agent:** BORDEN LADNER GERVAIS LLP

(54) **Titre : PROCÉDES DE SYNTHÈSE D'UN ANALOGUE DE PROSTACYCLINE**

(54) **Title: METHODS OF SYNTHESIZING A PROSTACYCLIN ANALOG**



IA

(57) **Abrégé/Abstract:**

The present invention provides processes for preparing a prostacyclin analogue of Formula (I) or a pharmaceutically acceptable salt thereof, wherein R¹⁰ is a linear or branched C₁₋₆ alkyl. The processes of the present invention comprise steps that generate

(57) Abrégé(suite)/Abstract(continued):

improved yields and fewer byproducts than traditional methods. The processes of the present invention employ reagents (e.g., the oxidizing reagent) that are less toxic than those used in the traditional methods (e.g., oxalyl chloride). Many of the processes of the present invention generate intermediates with improved e.e. and chemical purity; thereby eliminating the need of additional chromatography steps. And, the processes of the present invention are scalable to generate commercial quantities of the final compound.