



(86) Date de dépôt PCT/PCT Filing Date: 2009/10/16  
(87) Date publication PCT/PCT Publication Date: 2011/04/21  
(85) Entrée phase nationale/National Entry: 2012/04/10  
(86) N° demande PCT/PCT Application No.: US 2009/061076  
(87) N° publication PCT/PCT Publication No.: 2011/046569

(51) Cl.Int./Int.Cl. *C07D 307/87* (2006.01),  
*A61K 31/34* (2006.01), *A61P 9/02* (2006.01),  
*C07D 307/88* (2006.01), *C07D 307/885* (2006.01)  
(71) Demandeur/Applicant:  
CAYMAN CHEMICAL COMPANY, INCORPORATED,  
US  
(72) Inventeurs/Inventors:  
CHAMBOURNIER, GILLES, US;  
KORNILOV, ANDRIY, US;  
MAHMOUD, HUSSEIN M., US;  
VESELY, IVAN, CZ;  
BARRETT, STEPHEN DOUGLAS, US  
(74) Agent: GOWLING LAFLEUR HENDERSON LLP

(54) Titre : PROCEDE DE PREPARATION DE PROSTAGLANDINES DE LA SERIE F

(54) Title: PROCESS FOR THE PREPARATION OF F-SERIES PROSTAGLANDINS

(57) Abrégé/Abstract:

A process for the synthesis and purification of F-series prostaglandin compounds and synthetic intermediates used to prepare them. The synthetic intermediates are solid and may be purified by precipitation and therefore may form the representative F-series prostaglandin compounds such as latanoprost, bimatoprost, fluprostenol, cloprostenol, and substituted analogs therefrom in highly pure forms.