

# PRODUCT INFORMATION

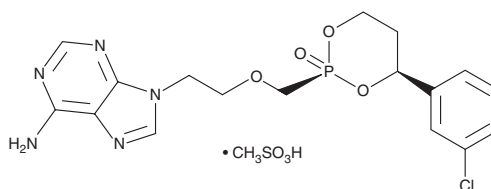


## Pradefovir (mesylate)

Item No. 45047

**CAS Registry No.:** 625095-61-6  
**Formal Name:** 9-[2-[[[(2R,4S)-4-(3-chlorophenyl)-2-oxido-1,3,2-dioxaphosphorinan-2-yl]methoxy]ethyl]-9H-purin-6-amine, methanesulfonate

**Synonym:** MB06866  
**MF:** C<sub>17</sub>H<sub>19</sub>ClN<sub>5</sub>O<sub>4</sub>P • CH<sub>3</sub>SO<sub>3</sub>H  
**FW:** 519.9  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Pradefovir (mesylate) is supplied as a solid. A stock solution may be made by dissolving the pradefovir (mesylate) in the solvent of choice, which should be purged with an inert gas. Pradefovir (mesylate) is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

### Description

Pradefovir is a liver-activated prodrug form of the antiviral nucleoside analog adefovir (Item No. 18650).<sup>1</sup> It is converted to adefovir diphosphate in the liver in rats when administered at a dose of 30 mg/kg.

### Reference

1. Reddy, K.R., Matelich, M.C., Ugarkar, B.G., *et al.* Pradefovir: A prodrug that targets adefovir to the liver for the treatment of hepatitis B. *J. Med. Chem.* **51**(3), 666-676 (2008).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/05/2026

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM