

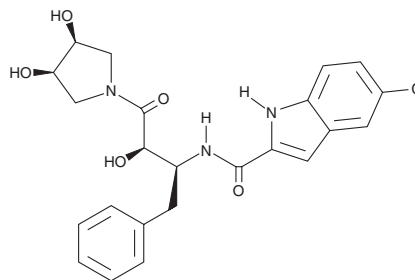
# PRODUCT INFORMATION



## Ingliforib

Item No. 38419

**CAS Registry No.:** 186392-65-4  
**Formal Name:** 5-chloro-N-[(1S,2R)-3-[(3R,4S)-3,4-dihydroxy-1-pyrrolidinyl]-2-hydroxy-3-oxo-1-(phenylmethyl)propyl]-1H-indole-2-carboxamide  
**Synonym:** CP 368,296  
**MF:** C<sub>23</sub>H<sub>24</sub>ClN<sub>3</sub>O<sub>5</sub>  
**FW:** 457.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 298 nm  
**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

Ingliforib is supplied as a solid. A stock solution may be made by dissolving the ingliforib in the solvent of choice, which should be purged with an inert gas. Ingliforib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ingliforib in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

### Description

Ingliforib is an inhibitor of the liver, muscle, and brain forms of glycogen phosphorylase (IC<sub>50</sub>s = 52, 352, and 150 nM, respectively).<sup>1</sup> It decreases myocardial glycogen phosphorylase activity and infarct size as a percentage of the area at risk in a rabbit model of ischemia-reperfusion injury induced by coronary artery occlusion.

### Reference

1. Tracey, W.R., Treadway, J.L., Magee, W.P., *et al.* Cardioprotective effects of ingliforib, a novel glycogen phosphorylase inhibitor. *Am. J. Physiol. Heart Circ. Physiol.* **286(3)**, H1177-H1184 (2004).

#### 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, 08/28/2023

#### 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