

# PRODUCT INFORMATION



**PF-06928215**

Item No. 32515

**CAS Registry No.:** 2378173-15-8  
**Formal Name:** (1R,2S)-2-[[[(7-hydroxy-5-phenylpyrazolo[1,5-a]pyrimidin-3-yl)carbonyl]amino]-cyclohexanecarboxylic acid

**MF:** C<sub>20</sub>H<sub>20</sub>N<sub>4</sub>O<sub>4</sub>

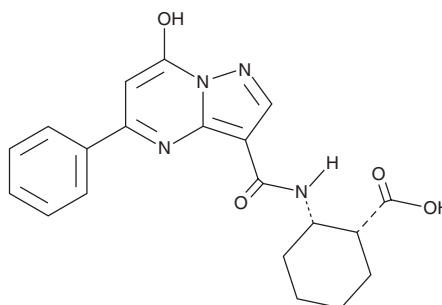
**FW:** 380.4

**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

PF-06928215 is supplied as a solid. A stock solution may be made by dissolving the PF-06928215 in the solvent of choice, which should be purged with an inert gas. PF-06928215 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

## Description

PF-06928215 is an inhibitor of cyclic GMP-AMP synthase (cGAS; IC<sub>50</sub> = 4.9 μM in a fluorescence polarization assay).<sup>1</sup>

## Reference

1. Hall, J., Brault, A., Vincent, F., *et al.* Discovery of PF-06928215 as a high affinity inhibitor of cGAS enabled by novel fluorescence polarization assay. *PLoS One* **12(9)**, e0184843 (2017).

### 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, 11/02/2022

## 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