

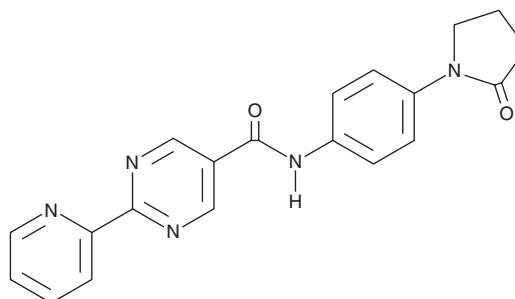
PRODUCT INFORMATION



Prostaglandin D Synthase (hematopoietic-type) Inhibitor F092

Item No. 28721

CAS Registry No.: 2250261-59-5
Formal Name: N-[4-(2-oxo-1-pyrrolidinyl)phenyl]-2-(2-pyridinyl)-5-pyrimidinecarboxamide
MF: C₂₀H₁₇N₅O₂
FW: 359.4
Purity: ≥98%
UV/Vis.: λ_{max}: 256, 285 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Prostaglandin D synthase (hematopoietic-type) inhibitor F092 is supplied as a crystalline solid. A stock solution may be made by dissolving the prostaglandin D synthase (hematopoietic-type) inhibitor F092 in the solvent of choice, which should be purged with an inert gas. Prostaglandin D synthase (hematopoietic-type) inhibitor F092 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of prostaglandin D synthase (hematopoietic-type) inhibitor F092 in these solvents is approximately 5 and 3 mg/ml, respectively. Prostaglandin D synthase (hematopoietic-type) inhibitor F092 is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of prostaglandin D synthase (hematopoietic-type) inhibitor F092 can be prepared by directly dissolving the crystalline solid in aqueous buffers. Prostaglandin D synthase (hematopoietic-type) inhibitor F092 is slightly soluble in PBS, pH 7.2. We do not recommend storing the aqueous solution for more than one day.

Description

Prostaglandin D synthase (hematopoietic-type) inhibitor F092 is an inhibitor of hematopoietic-type prostaglandin D synthase (H-PGDS; K_D = 0.14 nM).¹

Reference

1. Takaya, D., Inaka, K., Omura, A., *et al.* Characterization of crystal water molecules in a high-affinity inhibitor and hematopoietic prostaglandin D synthase complex by interaction energy studies. *Bioorg. Med. Chem.* **26(16)**, 4726-4734 (2018).

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, 04/01/2020

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