

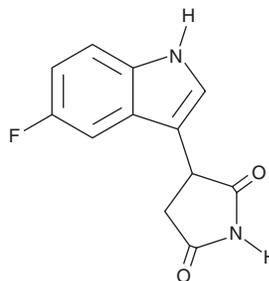
PRODUCT INFORMATION



EOS200271

Item No. 25003

CAS Registry No.: 198474-05-4
Formal Name: 3-(5-fluoro-1H-indol-3-yl)-2,5-pyrrolidinedione
Synonym: PF-06840003
MF: C₁₂H₉FN₂O₂
FW: 232.2
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 287 nm
Supplied as: A crystalline 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

EOS200271 is supplied as a crystalline solid. A stock solution may be made by dissolving the EOS200271 in the solvent of choice, which should be purged with an inert gas. EOS200271 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of EOS200271 in these solvents is approximately 30 mg/ml.

EOS200271 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, EOS200271 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. EOS200271 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

EOS200271 is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO-1; IC₅₀ = 0.15 μM for the human enzyme).¹ It is selective for IDO-1 over IDO-2 up to a concentration of 200 μM for the human recombinant enzymes. EOS200271 inhibits IDO-1-mediated production of kynurenine induced by IFN-γ in HeLa cells (IC₅₀ = 1.8 μM) and by LPS and IFN-γ in THP-1 cells and human-derived whole blood *ex vivo* (IC₅₀s = 1.7 and 4.7 μM, respectively).

Reference

1. Crosignani, S., Bingham, P.M., Bottemanne, P., *et al.* Discovery of a novel and selective indoleamine 2,3-dioxygenase (IDO-1) inhibitor 3-(5-fluoro-1H-indol-3-yl)pyrrolidine-2,5-dione (EOS200271/ PF-06840003) and its characterization as a potential clinical candidate. *J. Med. Chem.* 60(23), 9617-9629 (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, 12/19/2022

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