

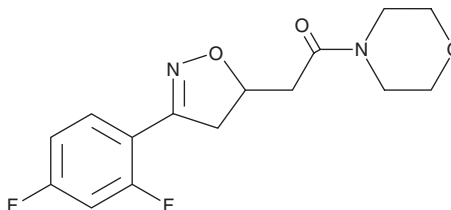
PRODUCT INFORMATION



CPSI-1306

Item No. 29905

CAS Registry No.: 1309793-47-2
Formal Name: 2-[3-(2,4-difluorophenyl)-4,5-dihydro-5-isoxazolyl]-1-(4-morpholinyl)-ethanone
MF: C₁₅H₁₆F₂N₂O₃
FW: 310.3
Purity: ≥95%
UV/Vis.: λ_{max}: 258 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

CPSI-1306 is supplied as a crystalline solid. A stock solution may be made by dissolving the CPSI-1306 in the solvent of choice, which should be purged with an inert gas. CPSI-1306 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CPSI-1306 in these solvents is approximately 30 mg/ml.

CPSI-1306 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPSI-1306 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CPSI-1306 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CPSI-1306 is an inhibitor of macrophage inhibitory factor (MIF).^{1,2} *In vivo*, CPSI-1306 (20 mg/kg per day) decreases skin thickness and myeloperoxidase (MPO) activity and induces keratinocyte apoptosis, as well as reduces papilloma formation and progression to micro-invasive squamous cell carcinoma (SCC) in a mouse model of UVB-induced SCC.¹ It lowers blood glucose levels and serum levels of IL-6 and TNF-α in a mouse model of non-insulin-dependent diabetes mellitus (NIDDM) induced by streptozotocin (Item No. 13104).²

References

1. Nagarajan, P., Tober, K.L., Riggenbach, J.A., *et al.* MIF antagonist (CPSI-1306) protects against UVB-induced squamous cell carcinoma. *Mol. Cancer Res.* **12(9)**, 1292-1302 (2014).
2. Sanchez-Zamora, Y., Terrazas, L.I., Vilches-Flores, A., *et al.* Macrophage migration inhibitory factor is a therapeutic target in treatment of non-insulin-dependent diabetes mellitus. *FASEB J.* **24(7)**, 2583-2590 (2010).

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, 11/16/2022

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