

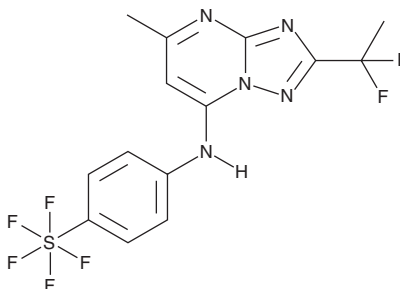
PRODUCT INFORMATION



DSM 265

Item No. 30680

CAS Registry No.: 1282041-94-4
Formal Name: (OC-6-21)-[4-[[2-(1,1-difluoroethyl)-5-methyl[1,2,4]triazolo[1,5-a]pyrimidin-7-yl]amino]phenyl]pentafluoro-sulfur
MF: $C_{14}H_{12}F_7N_5S$
FW: 415.3
Purity: $\geq 98\%$
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

DSM 265 is supplied as a solid. A stock solution may be made by dissolving the DSM 265 in the solvent of choice, which should be purged with an inert gas. DSM 265 is soluble in DMSO.

Description

DSM 265 is an antimalarial agent and inhibitor of *P. falciparum* and *P. vivax* dihydroorotate dehydrogenase (DHODH; $IC_{50}s = 8.9$ and 27 nM, respectively).¹ It is selective for *P. falciparum* and *P. vivax* DHODH over human, rabbit, pig, and monkey DHODH ($IC_{50}s = >41$ $\mu\text{g/ml}$) but does inhibit *P. cynomolgi*, dog, mouse, and rat DHODH ($IC_{50}s = 0.005$, 11 , 1.1 , and 0.82 $\mu\text{g/ml}$, respectively).² DSM 265 inhibits the growth of the *P. falciparum* strain 3D7 with an EC_{50} value of 4.3 nM.¹ It also reduces blood parasitemia in mice ($ED_{90} = 1.5$ mg/kg twice per day).²

References

1. Kokkonda, S., Deng, X., White, K.L., et al. Tetrahydro-2-naphthyl and 2-indanyl triazolopyrimidines targeting *Plasmodium falciparum* dihydroorotate dehydrogenase display potent and selective antimalarial activity. *J. Med. Chem.* **59**(11), 5416-5431 (2016).
2. Phillips, M.A., Lotharius, J., Marsh, K., et al. A long-duration dihydroorotate dehydrogenase inhibitor (DSM265) for prevention and treatment of malaria. *Sci. Transl. Med.* **7**(296), 296ra111 (2015).

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, 11/03/2022

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