

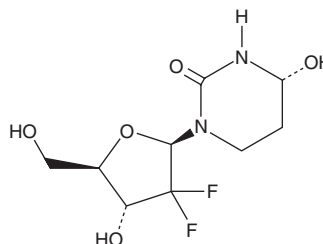
PRODUCT INFORMATION



Cedazuridine

Item No. 38767

CAS Registry No.: 1141397-80-9
Formal Name: (4R)-2'-deoxy-2',2'-difluoro-3,4,5,6-tetrahydro-uridine
Synonym: E7727
MF: C₉H₁₄F₂N₂O₅
FW: 268.2
Purity: ≥95%
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

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

Description

Cedazuridine is an inhibitor of cytidine deaminase ($IC_{50} = 0.4 \mu M$).¹ It increases plasma levels of decitabine via inhibition of its metabolism by cytidine deaminase in rhesus monkeys when administered at doses of 1 and 10 mg/kg. Cedazuridine (3 mg/kg) induces decreases in leukemic cell expansion in combination with orally administered 5-azacytidine (Item No. 11164), which is inactive as a single agent, in the bone marrow and spleen in a MOLM-13 acute myeloid leukemia (AML) mouse xenograft model.² Formulations containing cedazuridine in combination with decitabine have been used in the treatment of myelodysplastic syndrome (MDS).

References

1. Ferraris, D., Duvall, B., Delahanty, G., *et al.* Design, synthesis, and pharmacological evaluation of fluorinated tetrahydrouridine derivatives as inhibitors of cytidine deaminase. *J. Med. Chem.* **57**(6), 2582-2588 (2014).
2. Ramsey, H.E., Oganessian, A., Gorska, A.E., *et al.* Oral azacitidine and cedazuridine approximate parenteral azacitidine efficacy in murine model. *Target Oncol.* **15**(2), 231-240 (2020).

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, 09/18/2024

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