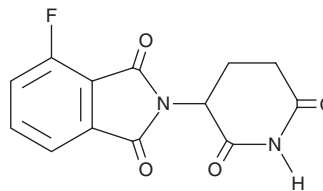


PRODUCT INFORMATION



TC-E3 5032
Item No. 26147

CAS Registry No.: 835616-60-9
Formal Name: 2-(2,6-dioxo-3-piperidiny)-4-fluoro-1H-isoindole-1,3(2H)-dione
MF: C₁₃H₉FN₂O₄
FW: 276.2
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 232, 240, 301 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

TC-E3 5032 is supplied as a crystalline solid. A stock solution may be made by dissolving the TC-E3 5032 in the solvent of choice, which should be purged with an inert gas. TC-E3 5032 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of TC-E3 5032 in these solvents is approximately 10 and 12 mg/ml, respectively.

TC-E3 5032 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TC-E3 5032 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TC-E3 5032 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TC-E3 5032 is a precursor in the synthesis of homodimeric proteolysis-targeting chimera technologies (PROTACs).¹ It has been used to attach various linkers at position 4 to produce alkylated pomalidomide derivatives as intermediates in the synthesis of homodimeric PROTACs. It has also been used as an intermediate in the synthesis of HDAC6 and anaplastic lymphoma kinase (ALK) degraders.^{2,3}

References

1. Steinebach, C., Lindner, S., Udeshi, N.D., *et al.* Homo-PROTACs for the chemical knockdown of cereblon. *ACS Chem. Biol.* **13**(9), 2771-2782 (2018).
2. Yang, K., Song, Y., Xie, H., *et al.* Development of the first small molecule histone deacetylase 6 (HDAC6) degraders. *Bioorg. Med. Chem. Lett.* **28**(14), 2493-2497 (2018).
3. Powell, C.E., Gao, Y., Tan, L., *et al.* Chemically induced degradation of anaplastic lymphoma kinase (ALK). *J. Med. Chem.* **61**(9), 4249-4255 (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, 11/28/2022

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