

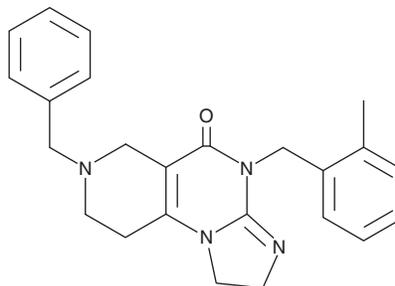
PRODUCT INFORMATION



TIC10

Item No. 16109

CAS Registry No.: 1616632-77-9
Formal Name: 2,4,6,7,8,9-hexahydro-4-[(2-methylphenyl)methyl]-7-(phenylmethyl)-imidazo[1,2-a]pyrido[3,4-e]pyrimidin-5(1H)-one
Synonym: ONC201
MF: C₂₄H₂₆N₄O
FW: 386.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 311 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

TIC10 is supplied as a crystalline solid. A stock solution may be made by dissolving the TIC10 in the solvent of choice, which should be purged with an inert gas. TIC10 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TIC10 in these solvents is approximately 10, 1, and 20 mg/ml, respectively.

TIC10 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TIC10 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TIC10 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

TIC10 induces the expression of TNF-related apoptosis-inducing ligand (TRAIL) at 1-5 μM in a p53-independent manner.¹ It is orally active, stable, and crosses the blood-brain barrier.¹ The induction of TRAIL results from dual inhibition of Akt and ERK1/2 combined with nuclear translocation of the transcription factor FOXO3a.¹ TIC10 suppresses the growth of orthotopic human glioblastoma multiforme tumors in mice.¹

Reference

1. Allen, J.E., Krigsfeld, G., Mayes, P.A., *et al.* Dual inactivation of Akt and ERK by TIC10 signals Foxo3a nuclear translocation, TRAIL gene induction, and potent antitumor effects. *Sci. Transl. Med.* **5(171)**, 171ra17 (2013).

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, 12/20/2022

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