

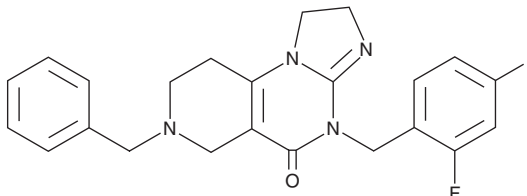
# PRODUCT INFORMATION



## ONC206

Item No. 36777

**CAS Registry No.:** 1638178-87-6  
**Formal Name:** 4-[(2,4-difluorophenyl)methyl]-2,4,6,7,8,9-hexahydro-7-(phenylmethyl)-imidazo[1,2-a]pyrido[3,4-e]pyrimidin-5(1H)-one  
**MF:** C<sub>23</sub>H<sub>22</sub>F<sub>2</sub>N<sub>4</sub>O  
**FW:** 408.4  
**Purity:** ≥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

ONC206 is supplied as a solid. A stock solution may be made by dissolving the ONC206 in the solvent of choice, which should be purged with an inert gas. ONC206 is soluble in DMSO and methanol.

### Description

ONC206 is a derivative of the TRAIL-inducing compound TIC10 (Item No. 16109) with anticancer activity.<sup>1</sup> It induces surface expression of TNF-related apoptosis-inducing ligand (TRAIL), apoptosis, and cell cycle arrest at the G<sub>1</sub> phase in HCT116 cells when used at a concentration of 0.05 μM. ONC206 is cytotoxic to six patient-derived diffuse intrinsic pontine glioma (DIPG) cell lines (IC<sub>50</sub>s = 0.05-0.173 μM).<sup>2</sup>

### References

1. Wagner, J., Kline, C.L., Ralff, M.D., *et al.* Preclinical evaluation of the imipridone family, analogs of clinical stage anti-cancer small molecule ONC201, reveals potent anti-cancer effects of ONC212. *Cell Cycle* **16(19)**, 1790-1799 (2017).
2. Borsuk, R., Zhou, L., Chang, W.-I., *et al.* Potent preclinical sensitivity to imipridone-based combination therapies in oncohistone H3K27M-mutant diffuse intrinsic pontine glioma is associated with induction of the integrated stress response, TRAIL death receptor DR5, reduced ClpX and apoptosis. *Am. J. Cancer Res.* **11(9)**, 4607-4623 (2021).

#### 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, 02/21/2024

#### 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