

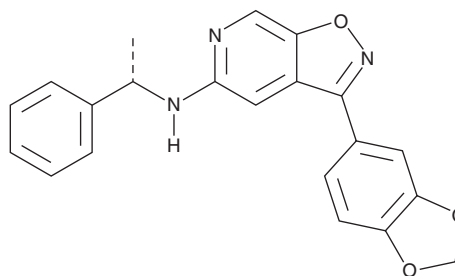
# PRODUCT INFORMATION



## TC-S 7005

Item No. 45232

**CAS Registry No.:** 1082739-92-1  
**Formal Name:** 3-(1,3-benzodioxol-5-yl)-N-[(1S)-1-phenylethyl]-isoxazolo[5,4-c]pyridin-5-amine  
**MF:** C<sub>21</sub>H<sub>17</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 359.4  
**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

TC-S 7005 is supplied as a solid. A stock solution may be made by dissolving the TC-S 7005 in the solvent of choice, which should be purged with an inert gas. TC-S 7005 is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

TC-S 7005 is an inhibitor of polo-like kinase 2 (Plk2; IC<sub>50</sub> = 4 nM).<sup>1</sup> It is selective for Plk2 over Plk1 and Plk3 in a cell-free assay (IC<sub>50</sub>s = 214 and 24 nM, respectively). It induces mitotic arrest and cell death in HCT116 colorectal cancer cells. TC-S 7005 reduces IL-1β, IL-6, inducible nitric-oxide synthase (iNOS), and COX-2 levels in LPS-stimulated BV-2 microglial cells when used at a concentration of 8 μM and decreases NF-κB transcriptional activation in the same cells.<sup>2</sup>

### References

- Hanan, E.J., Fucini, R.V., Romanowski, M.J., *et al.* Design and synthesis of 2-amino-isoxazopyridines as Polo-like kinase inhibitors. *Bioorg. Med. Chem. Lett.* **18**(19), 5186-5189 (2008).
- Cheng, J., Wu, L., Chen, X., *et al.* Polo-like kinase 2 promotes microglial activation via regulation of the HSP90α/IKKβ pathway. *Cell Rep.* **43**(10), 114827 (2024).

#### 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, 02/27/2026

#### 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