

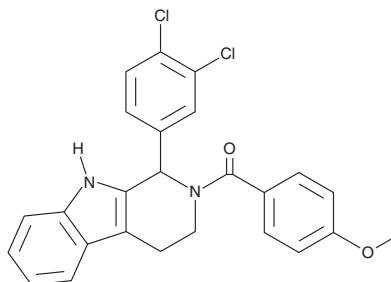
# PRODUCT INFORMATION



## CAY10719

Item No. 20469

**CAS Registry No.:** 1942919-63-2  
**Formal Name:** [1-(3,4-dichlorophenyl)-1,3,4,9-tetrahydro-2H-pyrido[3,4-b]indol-2-yl](4-methoxyphenyl)-methanone  
**MF:** C<sub>25</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 451.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 225 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

CAY10719 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10719 in the solvent of choice. CAY10719 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CAY10719 in these solvents is approximately 2 and 20 mg/ml, respectively.

CAY10719 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10719 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10719 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CAY10719 is a selective inhibitor of the breast cancer resistance protein ABCG2 (IC<sub>50</sub> = 0.23 μM) with little activity at ABCG1.<sup>1</sup> It has been shown to reverse the ABCG2-mediated resistance toward SN 38 (Item No. 15632) and to inhibit ATPase activity.<sup>1</sup>

### Reference

1. Spindler, A., Stefan, K. and Wiese, M. Synthesis and investigation of tetrahydro-β-carboline derivatives as inhibitors of the breast cancer resistance protein (ABCG2). *J. Med. Chem.* **59(13)**, 6121-6135 (2016).

#### 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, 08/08/2022

#### 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