

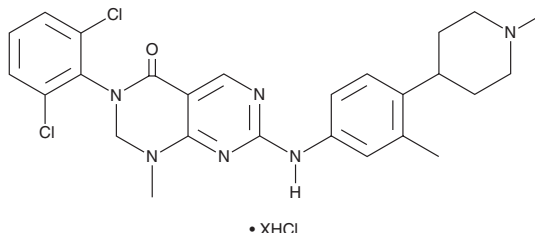
# PRODUCT INFORMATION



## Debio 0123 (hydrochloride)

Item No. 39674

**Formal Name:** 3-(2,6-dichlorophenyl)-2,3-dihydro-1-methyl-7-[[3-methyl-4-(1-methyl-4-piperidiny)phenyl]amino]-pyrimido[4,5-d]pyrimidin-4(1H)-one, hydrochloride  
**Synonyms:** WEE1-IN-5, WEE1 Inhibitor 5  
**MF:** C<sub>26</sub>H<sub>28</sub>Cl<sub>2</sub>N<sub>6</sub>O • XHCl  
**FW:** 511.4  
**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

Debio 0123 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the debio 0123 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Debio 0123 (hydrochloride) is soluble in methanol.

### Description

Debio 0123 is an inhibitor of the checkpoint kinase WEE1 (IC<sub>50</sub> = 0.8 nM).<sup>1</sup> *In vivo*, debio 0123 (30 mg/kg per day), in combination with the inhibitor of topoisomerase II etoposide (Item No. 12092), reduces tumor volume in a patient-derived xenograft (PDX) mouse model of stomach cancer.<sup>2</sup>

### References

1. Hewitt, P., Burkamp, F., Wilkinson, A., *et al.* Pyrimidopyrimidones useful as WEE-1 kinase inhibitors. *Debiopharm International S.A. WO 2023/170065 A1* (2018).
2. Rodrigo Imedio, E., Piggott, L., Bellon, A., *et al.* Methods of treating small cell lung cancer. *Almac Discovery Limited. WO 2018/162932 A1* (2023).

#### 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, 08/22/2024

#### 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](http://WWW.CAYMANCHEM.COM)