

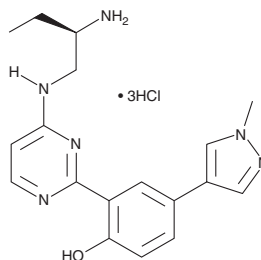
# PRODUCT INFORMATION



## CRT0066101 (hydrochloride)

Item No. 15337

**CAS Registry No.:** 1781742-22-0  
**Formal Name:** 2-[4-[[[(2R)-2-aminobutyl]amino]-2-pyrimidinyl]-4-(1-methyl-1H-pyrazol-4-yl)-phenol, trihydrochloride  
**MF:** C<sub>18</sub>H<sub>22</sub>N<sub>6</sub>O • 3HCl  
**FW:** 447.8  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 250 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

CRT0066101 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CRT0066101 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CRT0066101 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CRT0066101 (hydrochloride) in these solvents is approximately 3 and 0.1 mg/ml, respectively.

CRT0066101 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CRT0066101 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CRT0066101 (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Protein kinase D (PKD) is a serine/threonine protein kinase that is activated by diacylglycerol, commonly downstream of PKC signaling. The three human PKD isoforms target a variety of proteins to alter cell proliferation, survival, invasion, and protein transport. CRT0066101 is an inhibitor of all three PKD isoforms (IC<sub>50</sub>s = 1, 2.5, and 2 nM for PKD1, PKD2, and PKD3, respectively).<sup>1</sup> It exhibits selectivity for PKD against a panel of >90 protein kinases, including PKCα, PKBα, MEK, ERK, c-Raf, c-Src, and c-Abl.<sup>1</sup> It can block cell proliferation, induce apoptosis, and reduce the viability of pancreatic cancer cells both *in vitro* and *in vivo*.<sup>1</sup>

### Reference

1. Harikumar, K.B., Kunnumakkara, A.B., Ochi, N., *et al.* A novel small-molecule inhibitor of protein kinase D blocks pancreatic cancer growth *in vitro* and *in vivo*. *Mol. Cancer Ther.* **9**(5), 1136-1146 (2010).

#### 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, 12/21/2022

#### 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