

PRODUCT INFORMATION



CAY10740 (hydrochloride)

Item No. 28320

Formal Name: ((3S,4R)-3-amino-4-hydroxypiperidin-1-yl) (2-(1-benzyl-1H-indol-2-yl)-7-methoxy-1-methyl-1H-benzo[d]imidazol-5-yl)methanone, monohydrochloride

MF: C₃₀H₃₁N₅O₃ • HCl

FW: 546.1

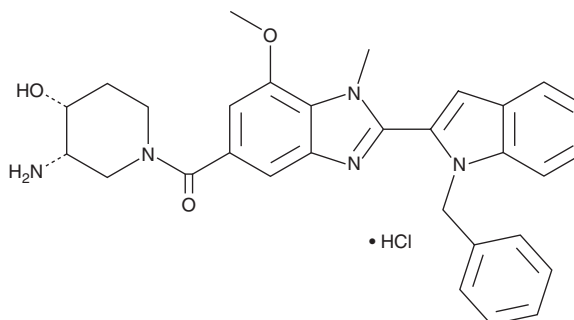
Purity: ≥98%

UV/Vis.: λ_{max}: 244, 304 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

CAY10740 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10740 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CAY10740 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10740 (hydrochloride) in these solvents is approximately 0.3, 5, and 10 mg/ml, respectively.

CAY10740 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10740 (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10740 (hydrochloride) has a solubility of approximately 0.3 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

CAY10740 is an analog of the protein arginine deiminase 4 (PAD4) inhibitor GSK484 (Item No. 17488).¹ It decreases the viability of 4T1 cells by 85 and 98% when used at concentrations of 5 and 10 μM, respectively, which is more than five-fold as effective as GSK484.

Reference

1. Guo, Z., Shi, L., Wang, B., *et al.* Synthesis of reversible PAD4 inhibitors via copper-catalyzed C–H arylation of benzimidazole. *Sci. China Chem.* **62**(6), 592-596 (2019).

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, 05/15/2020

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