

PRODUCT INFORMATION

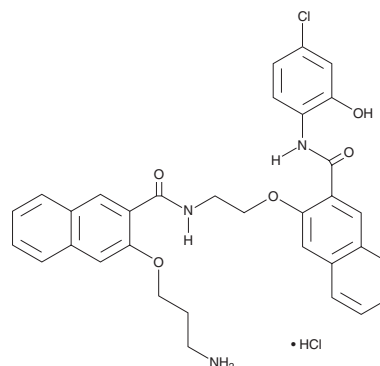


666-15

Item No. 30780

CAS Registry No.: 1433286-70-4
Formal Name: 3-(3-aminopropoxy)-N-[2-[[[3-[[[4-chloro-2-hydroxyphenyl]amino]carbonyl]-2-naphthalenyl]oxy]ethyl]-2-naphthalenecarboxamide, monohydrochloride

Synonym: Compound 3i
MF: C₃₃H₃₀ClN₃O₅ • HCl
FW: 620.5
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm
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

666-15 is supplied as a solid. A stock solution may be made by dissolving the 666-15 in the solvent of choice, which should be purged with an inert gas. 666-15 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 666-15 in these solvents is approximately 5 mg/ml.

Description

666-15 is an inhibitor of gene transcription mediated by cAMP-response-element-binding protein (CREB; IC₅₀ = 81 nM in a cell-based reporter assay).¹ It inhibits the growth of A549, MCF-7, MDA-MB-231, and MDA-MB-468 cells (GI₅₀s = 0.47, 0.31, 0.073, and 0.046 μM, respectively). 666-15 (10 mg/kg per day, five days per week) reduces tumor growth in an MDA-MB-468 mouse xenograft model.

Reference

1. Xie, F., Li, B.X., Kassenbrock, A., *et al.* Identification of a potent inhibitor of CREB-mediated gene transcription with efficacious in vivo anticancer activity. *J. Med. Chem.* **58**(12), 5075-5087 (2015).

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, 12/20/2022

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