

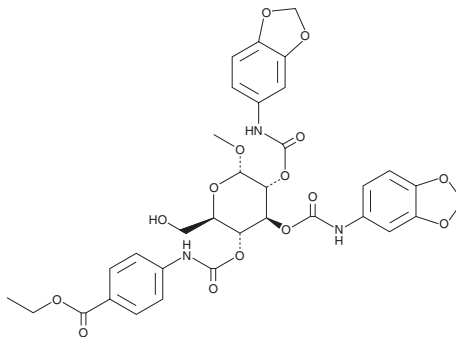
PRODUCT INFORMATION



BRD7552

Item No. 19143

CAS Registry No.: 1137359-47-7
Formal Name: α -D-glucopyranoside, methyl, 2,3-bis(N-1,3-benzodioxol-5-ylcarbamate) 4-[N-[4-(ethoxycarbonyl)phenyl]carbamate]
MF: C₃₃H₃₃N₃O₁₅
FW: 711.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 252 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

BRD7552 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD7552 in the solvent of choice, which should be purged with an inert gas. BRD7552 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD7552 in these solvents is approximately 10, 20, and 25 mg/ml, respectively.

BRD7552 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD7552 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD7552 has a solubility of approximately 0.2 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD7552 is a small molecule that induces expression of pancreatic and duodenal homeobox 1 (PDX1), a transcription factor involved in pancreas development and function.¹ BRD7552 dose-dependently increases PDX1 and insulin gene expression, as well as PDX1 protein levels, in PANC-1 cells. It also increases acetylation of histone H3 and trimethylation of H3K4 and H3K9.

Reference

1. Yuan, Y., Hartland, K., Boxkovic, Z., *et al.* A small-molecule inducer of PDX1 expression identified by high-throughput screening. *Chem. Biol.* **20**(12), 1513-1522 (2013).

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, 11/08/2022

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