

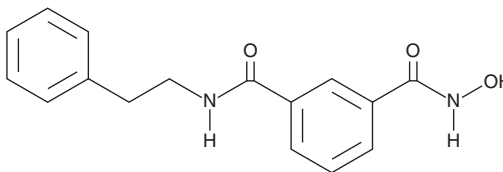
# PRODUCT INFORMATION



**BRD73954**

Item No. BRD73954

CAS Registry No.: 1440209-96-0  
Formal Name: N<sup>1</sup>-hydroxy-N<sup>3</sup>-(2-phenylethyl)-  
1,3-benzenedicarboxamide  
MF: C<sub>16</sub>H<sub>16</sub>N<sub>2</sub>O<sub>3</sub>  
FW: 284.3  
Purity: ≥95%  
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

BRD73954 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD73954 in the solvent of choice, which should be purged with an inert gas. BRD73954 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD73954 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

BRD73954 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD73954 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BRD73954 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

BRD73954 is a dual inhibitor of histone deacetylase 6 (HDAC6) and HDAC8 (IC<sub>50</sub>s = 36 and 120 nM, respectively).<sup>1</sup> It is selective for HDAC6 and -8 over HDAC1-5, -7, and -9 (IC<sub>50</sub>s = 12, 9, 23, >33, >33, 13, and >33 μM, respectively). BRD73954 (10 μM) increases acetylation of α-tubulin, a known HDAC6 substrate, but not histone H3, a substrate for HDAC1, -2, and -3, in HeLa cells.

## Reference

1. Olson, D.E., Wagner, F.F., Kaya, T., *et al.* Discovery of the first histone deacetylase 6/8 dual inhibitors. *J. Med. Chem.* **56**(11), 4816-4820 (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, 11/30/2022

## 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