

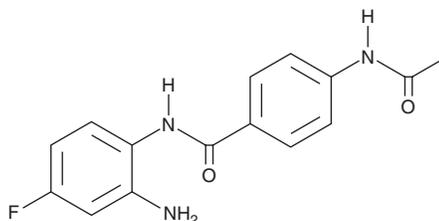
PRODUCT INFORMATION



BRD3308

Item No. 34875

CAS Registry No.: 1550053-02-5
Formal Name: 4-(acetlamino)-N-(2-amino-4-fluorophenyl)-benzamide
MF: C₁₅H₁₄FN₃O₂
FW: 287.3
Purity: ≥98%
UV/Vis.: λ_{max}: 243, 286, 303 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

BRD3308 is supplied as a solid. A stock solution may be made by dissolving the BRD3308 in the solvent of choice, which should be purged with an inert gas. BRD3308 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BRD3308 in these solvents is approximately 30 mg/ml.

Description

BRD3308 is an inhibitor of histone deacetylase 3 (HDAC3; K_i = 29 nM).¹ It is selective for HDAC3 over HDAC1 and HDAC2 (K_is = 6,300 and 5,100 nM, respectively). BRD3308 (10 μM) reduces cytokine-induced increases in apoptosis in rat INS-1E β cells. *In vivo*, BRD3308 (5 mg/kg) reduces plasma glucose levels in rats and increases plasma insulin levels in rats in a hyperglycemic clamp assay as a model of type 2 diabetes.²

References

1. Wagner, F.F., Lundh, M., Kaya, T., *et al.* An isochemogenic set of inhibitors To define the therapeutic potential of histone deacetylases in β-Cell protection. *ACS Chem. Biol.* **11**(2), 363-374 (2016).
2. Lundh, M., Galbo, T., Poulsen, S.S., *et al.* Histone deacetylase 3 inhibition improves glycaemia and insulin secretion in obese diabetic rats. *Diabetes Obes. Metab.* **17**(7), 703-707 (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, 10/17/2022

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