

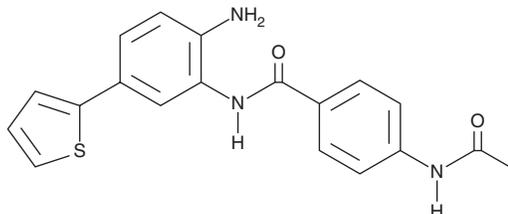
# PRODUCT INFORMATION



## BRD6929

Item No. 34149

**CAS Registry No.:** 849234-64-6  
**Formal Name:** 4-(acetilamino)-N-[2-amino-5-(2-thienyl)phenyl]-benzamide  
**Synonym:** JQ12  
**MF:** C<sub>19</sub>H<sub>17</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 351.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 282 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

BRD6929 is supplied as a solid. A stock solution may be made by dissolving the BRD6929 in the solvent of choice, which should be purged with an inert gas. BRD6929 is soluble in DMSO.

### Description

BRD6929 is a histone deacetylase 1 (HDAC1) and HDAC2 inhibitor (IC<sub>50</sub>s = 0.04 and 0.1 μM, respectively, for the recombinant human enzymes).<sup>1</sup> It is selective for HDAC1 and HDAC2 over HDAC3 and HDAC8 (IC<sub>50</sub>s = >20 μM for both). BRD6929 inhibits proliferation of HCT116 human colon cancer cells but not non-cancerous human mammary epithelial cells (HMECs; IC<sub>50</sub>s = 4 and >50 μM, respectively).

### Reference

1. Moradei, O.M., Mallais, T.C., Frechette, S., *et al.* Novel aminophenyl benzamide-type histone deacetylase inhibitors with enhanced potency and selectivity. *J. Med. Chem.* **50**(23), 5543-5546 (2007).

#### 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, 12/01/2022

#### 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