

# PRODUCT INFORMATION



**J27644**

Item No. 39469

**Formal Name:** 4-((8-chloro-2,4-dioxo-3,4-dihydroquinazolin-1(2H)-yl)methyl)-N-hydroxybenzamide

**MF:** C<sub>16</sub>H<sub>12</sub>ClN<sub>3</sub>O<sub>4</sub>

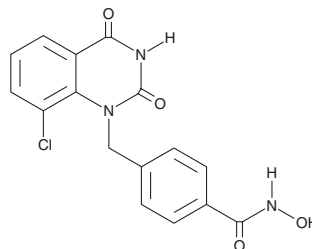
**FW:** 345.7

**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥3 years



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

## Laboratory Procedures

J27644 is supplied as a solid. A stock solution may be made by dissolving the J27644 in the solvent of choice, which should be purged with an inert gas. J27644 is soluble in DMSO and methanol.

## Description

J27644 is an inhibitor of histone deacetylase 6 (HDAC6; IC<sub>50</sub> = 10 nM).<sup>1</sup> It is selective for HDAC6 over HDAC1-5 (IC<sub>50</sub>s = 211-516 nM) and HDAC7, -8, -9, and -11 (IC<sub>50</sub>s = 137-812 nM). J27644 (10 μM) reduces the level of proteins involved in fibrosis, including collagen 1A1, α-smooth muscle actin (α-SMA), and fibronectin, in human pulmonary fibroblasts *in vitro*.

## Reference

1. Yu, W.-C., Yeh, T.-Y., Ye, C.-H., *et al.* Discovery of HDAC6, HDAC8, and 6/8 inhibitors and development of cell-based drug screening models for the treatment of TGF-β-induced idiopathic pulmonary fibrosis. *J. Med. Chem.* **66**(15), 10528-10557 (2023).

### 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, 10/26/2023

## 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