

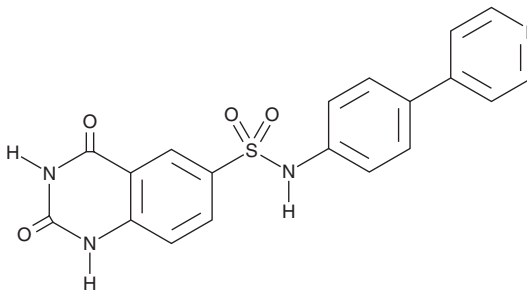
# PRODUCT INFORMATION



**ARN19874**

Item No. 24283

**CAS Registry No.:** 2190502-57-7  
**Formal Name:** 1,2,3,4-tetrahydro-2,4-dioxo-N-[4-(4-pyridinyl)phenyl]-6-quinazolinesulfonamide  
**MF:** C<sub>19</sub>H<sub>14</sub>N<sub>4</sub>O<sub>4</sub>S  
**FW:** 394.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220, 264 nm  
**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

ARN19874 is supplied as a crystalline solid. A stock solution may be made by dissolving the ARN19874 in the solvent of choice, which should be purged with an inert gas. ARN19874 is soluble in the organic solvent DMSO, with warming and sonication, at a concentration of approximately 1 mg/ml.

## Description

ARN19874 is an inhibitor of N-acyl phosphatidylethanolamine phospholipase D (NAPE-PLD; IC<sub>50</sub> = 33.7 μM).<sup>1</sup> It is selective for NAPE-PLD over carbonic anhydrase II, neutral endopeptidase, and angiotensin-converting enzyme (ACE) up to 50 μM. ARN19874 (50 μM) increases the levels of certain NAPE substrates and decreases the endogenous levels of stearoyl ethanolamide (Item No. 90245), but not oleoyl ethanolamide (Item No. 90265) or palmitoyl ethanolamide (Item No. 90350), in HEK293 cells.

## Reference

1. Castellani, B., Diamanti, E., Pizzirani, D., *et al.* Synthesis and characterization of the first inhibitor of N-acylphosphatidylethanolamine phospholipase D (NAPE-PLD). *Chem. Commun. (Camb.)* **53(95)**, 12814-12817 (2017).

### 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, 12/06/2022

## 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