

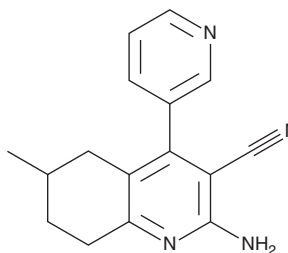
# PRODUCT INFORMATION



## BRD6989

Item No. 35358

**CAS Registry No.:** 642008-81-9  
**Formal Name:** 2-amino-5,6,7,8-tetrahydro-6-methyl-4-(3-pyridinyl)-3-quinolinecarbonitrile  
**MF:** C<sub>16</sub>H<sub>16</sub>N<sub>4</sub>  
**FW:** 264.3  
**Purity:** ≥98%  
**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

BRD6989 is supplied as a solid. A stock solution may be made by dissolving the BRD6989 in the solvent of choice, which should be purged with an inert gas. BRD6989 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

BRD6989 is an inhibitor of the Cdk8/cyclin C complex.<sup>1</sup> It selectively binds to the Cdk8/cyclin C complex (IC<sub>50</sub> = ~200 nM) over the Cdk1/cyclin B, Cdk2/cyclin A, Cdk5/p35, Cdk7/cyclin H, and Cdk9/cyclin T complexes at 1,000 nM. BRD6989 increases the production of IL-10 in primary mouse bone marrow-derived dendritic cells (BMDCs) stimulated with zymosan A (Item No. 21175) with an EC<sub>50</sub> value of approximately 1 μM and reduces zymosan A-induced increases in IL-6 levels in the same cells.

### Reference

1. Johannessen, L., Sundberg, T.B., O'Connell, D.J., *et al.* Small-molecule studies identify CDK8 as a regulator of IL-10 in myeloid cells. *Nat. Chem. Biol.* **13**(10), 1102-1108 (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, 12/12/2022

#### 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