

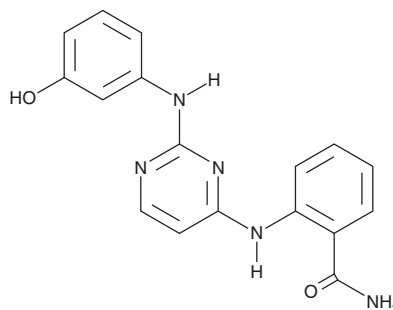
# PRODUCT INFORMATION



**DB07268**

Item No. 22257

**CAS Registry No.:** 929007-72-7  
**Formal Name:** 2-[[2-[(3-hydroxyphenyl)amino]-4-pyrimidinyl]amino]-benzamide  
**MF:** C<sub>17</sub>H<sub>15</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 321.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 275, 323 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

DB07268 is supplied as a crystalline solid. A stock solution may be made by dissolving the DB07268 in the solvent of choice. DB07268 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of DB07268 in these solvents is approximately 30 mg/ml.

DB07268 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DB07268 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DB07268 has a solubility of approximately 0.03 mg/ml in a 1:30 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

DB07268 is a potent inhibitor of JNK1 (IC<sub>50</sub> = 9 nM) that binds to the ATP site of JNK1.<sup>1</sup> It is selective for JNK1 over P38, ERK2, AKT1, CHK1, and PAK4 among others. DB07268 did not reduce phosphorylation of insulin resistance substrate 1 (IRS-1) in HEK293 cells.<sup>2</sup>

## References

1. Liu, M., Wang, S., Clampit, J.E., *et al.* Discovery of a new class of 4-anilinopyrimidines as potent c-Jun N-terminal kinase inhibitors: Synthesis and SAR studies. *Bioorg. Med. Chem. Lett.* **17(3)**, 668-672 (2007).
2. Simon-Szabó, S., Kokas, M., Greff, Z., *et al.* Novel compounds reducing IRS-1 serine phosphorylation for treatment of diabetes. *Bioorg. Med. Chem. Lett.* **26(2)**, 424-428 (2016).

### 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, 12/22/2022

## 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