

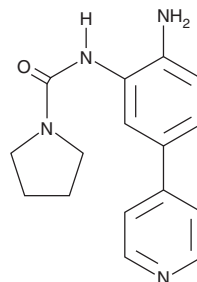
PRODUCT INFORMATION



BRD6688

Item No. 19836

CAS Registry No.: 1404562-17-9
Formal Name: N-[2-amino-5-(4-pyridinyl)phenyl]-1-pyrrolidinecarboxamide
MF: C₁₆H₁₈N₄O
FW: 282.3
Purity: ≥98%
UV/Vis.: λ_{max}: 315 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

BRD6688 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD6688 in the solvent of choice, which should be purged with an inert gas. BRD6688 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD6688 in ethanol is approximately 2 mg/ml and approximately 30 in DMSO and DMF.

BRD6688 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD6688 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD6688 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD6688 is an HDAC inhibitor with IC₅₀ values of 21 nM, 100 nM, and 11.48 μM for HDAC1, 2, and 3, respectively.¹ It possesses preferential binding kinetics with six-fold extended half-life on HDAC2 compared to HDAC1 (381 min versus 65 min).¹ BRD6688 crosses the blood brain barrier and has been shown to rescue the memory defects associated with p25 induced neurodegeneration in contextual fear conditioning in a CK-p25 mouse model of neurodegeneration.¹

Reference

1. Wagner, F. F., Zhang, Y.-L., Fass, D. M., *et al.* Kinetically selective inhibitors of histone deacetylase 2 (HDAC2) as cognition enhancers. *Chem. Sci.* **6**(1), 804-815 (2015).

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, 10/13/2022

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