

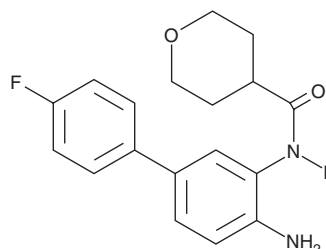
PRODUCT INFORMATION



BRD4884

Item No. 19834

CAS Registry No.: 1404559-91-6
Formal Name: N-(4-amino-4'-fluoro[1,1'-biphenyl]-3-yl)tetrahydro-2H-pyran-4-carboxamide
MF: C₁₈H₁₉FN₂O₂
FW: 314.4
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 276 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

BRD4884 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD4884 in the solvent of choice, which should be purged with an inert gas. BRD4884 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD4884 in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

BRD4884 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD4884 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BRD4884 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD4884 is an HDAC inhibitor with IC₅₀ values of 29 nM, 62 nM, and 1.09 μM for HDAC1, 2, and 3, respectively.¹ It possesses preferential binding kinetics with seven-fold longer half-life on HDAC2 compared to HDAC1 (143 min versus 20 min).¹ BRD4884 crosses the blood brain barrier and has been evaluated in CK-p25 mice, a 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, 12/09/2022

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