

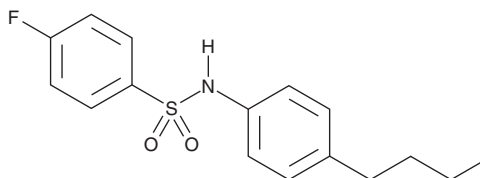
# PRODUCT INFORMATION



**DC260126**

Item No. 19388

**CAS Registry No.:** 346692-04-4  
**Formal Name:** N-(4-butylphenyl)-4-fluorobenzenesulfonamide  
**MF:** C<sub>16</sub>H<sub>18</sub>FNO<sub>2</sub>S  
**FW:** 307.4  
**Purity:** ≥98%  
**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

DC260126 is supplied as a crystalline solid. A stock solution may be made by dissolving the DC260126 in the solvent of choice. DC260126 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of DC260126 in these solvents is approximately 100, 20, and 30 mg/ml, respectively.

DC260126 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DC260126 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. DC260126 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

DC260126 is a free fatty acid receptor 1 (FFAR1/GPR40) antagonist that inhibits GPR40-mediated Ca<sup>2+</sup> elevations stimulated by linoleic, oleic, palmitoleic, and lauric acid (Item Nos. 90150, 90260, 10009871, and 10006626, respectively) with IC<sub>50</sub> values of 6.28, 5.96, 7.07, and 4.58 μM, respectively.<sup>1</sup> At 10 mg/kg, it has been shown to inhibit glucose-stimulated insulin secretion, to reduce blood insulin levels, to improve insulin sensitivity, and to decrease the rate of pancreatic β-cell apoptosis in obese diabetic *db/db* mice.<sup>2</sup>

## References

1. Hu, H., He, L., Gong, Z., *et al.* A novel class of antagonists for the FFAs receptor GPR40. *Biochem. Bioph. Res. Commun.* **390**(3), 557-563 (2009).
2. Sun, T., Wang, T., Zhou, Y., *et al.* DC260126: A small-molecule antagonist of GPR40 that protects against pancreatic β-cells dysfunction in *db/db* mice. *PLoS One* **8**(6), (2013).

### 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, 01/03/2023

## 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