# **PRODUCT** INFORMATION



HJC0350

Item No. 22949

CAS Registry No.:	885434-70-8	
Formal Name:	2,4-dimethyl-1-[(2,4,6-trimethylphenyl)sulfonyl]-1H-pyrrole	\
MF:	C <sub>15</sub> H <sub>19</sub> NO <sub>2</sub> S	
FW:	277.4	
Purity:	≥98%	
UV/Vis.:	λ <sub>max</sub> : 204 nm	N S
Supplied as:	A crystalline solid	
Storage:	-20°C	, 0.0
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

## Laboratory Procedures

HJC0350 is supplied as a crystalline solid. A stock solution may be made by dissolving the HJC0350 in the solvent of choice. HJC0350 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of HJC0350 in these solvents is approximately 10, 30, and 20 mg/ml, respectively.

HJC0350 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HJC0350 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HJC0350 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

HJC0350 is a selective inhibitor of exchange protein directly activated by cAMP 2 (Epac2) with an  $IC_{50}$  value of 300 nM for competition with 8-NBD-cAMP binding.<sup>1</sup> It inhibits Epac2 guanine nucleotide exchange activity but does not inhibit Epac1-mediated Rap1-GDP exchange or cAMP-mediated protein kinase A (PKA) activation in vitro, indicating it is selective for Epac2 at a concentration of 25  $\mu$ M. HJC0350 also blocks stimulation of Epac2, but not Epac1, by 007-AM in HEK293 cells.

## Reference

1. Chen, H., Tsalkova, T., Chepurny, O.G., et al. Identification and characterization of small molecules as potent and specific EPAC2 antagonists. J. Med. Chem. 56(3), 952-962 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFFTY 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

uyer 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, 10/04/2022

# 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