

PRODUCT INFORMATION

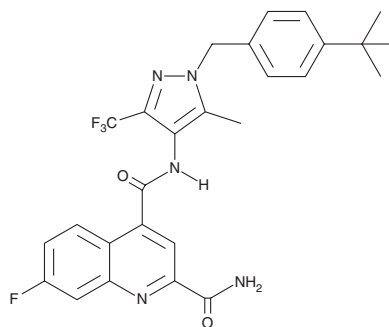


BAY-588

Item No. 19960

CAS Registry No.: 1799759-24-2
Formal Name: N⁴-[1-[4-(1,1-dimethylethyl)phenyl]methyl]-5-methyl-3-(trifluoromethyl)-1H-pyrazol-4-yl]-7-fluoro-2,4-quinolinedicarboxamide

MF: C₂₇H₂₅F₄N₅O₂
FW: 527.5
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 292 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

BAY-588 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY-588 in the solvent of choice, which should be purged with an inert gas. BAY-588 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BAY-588 in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

BAY-588 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY-588 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BAY-588 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

BAY-588 is a glucose transporter inhibitor and derivative of BAY-876 (Item No. 19961).¹ It selectively inhibits glucose transporter 4 (Glut4), Glut1, and Glut3 over Glut2 with IC₅₀ values of 0.5, 1.18, 5.47, and >10 μM, respectively, in CHO cells expressing human recombinant transporters.

Reference

1. Siebeneicher, H., Cleve, A., Rehwinkel, H., *et al.* Identification and optimization of the first highly selective GLUT1 inhibitor BAY-876. *ChemMedChem* **11**(20), 2261-2271 (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, 11/01/2022

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