

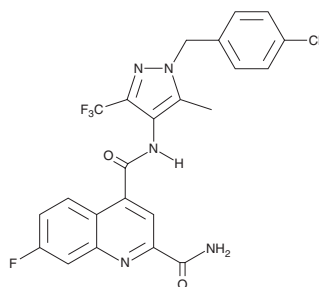
# PRODUCT INFORMATION



## BAY-876

Item No. 19961

**CAS Registry No.:** 1799753-84-6  
**Formal Name:** N<sup>4</sup>-[1-[(4-cyanophenyl)methyl]-5-methyl-3-(trifluoromethyl)-1H-pyrazol-4-yl]-7-fluoro-2,4-quinolinedicarboxamide  
**MF:** C<sub>24</sub>H<sub>16</sub>F<sub>4</sub>N<sub>6</sub>O<sub>2</sub>  
**FW:** 496.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234 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-876 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY-876 in the solvent of choice, which should be purged with an inert gas. BAY-876 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BAY-876 in these solvents is approximately 30 mg/ml.

BAY-876 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY-876 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BAY-876 has a solubility of approximately 0.33 mg/ml in a 1:2 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-876 is a potent and selective inhibitor of glucose transporter 1 (Glut1) with IC<sub>50</sub> values of 2, 10,800, 1,670, and 290 nM for Glut1, Glut2, Glut3, and Glut4, respectively, in CHO cells expressing human recombinant receptors.<sup>1</sup> It shows low metabolic clearance and has high permeability *in vitro*. *In vivo*, BAY-876 displays low plasma clearance and high oral bioavailability in rats and dogs.

### 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, 12/08/2022

#### 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