

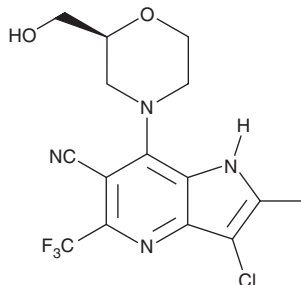
# PRODUCT INFORMATION



**PF-06869206**

Item No. 28158

**CAS Registry No.:** 2227425-05-8  
**Formal Name:** 3-chloro-7-[(2S)-2-(hydroxymethyl)-4-morpholinyl]-2-methyl-5-(trifluoromethyl)-1H-pyrrolo[3,2-b]pyridine-6-carbonitrile  
**MF:** C<sub>15</sub>H<sub>14</sub>ClF<sub>3</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 374.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 250, 282, 324 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

PF-06869206 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-06869206 in the solvent of choice, which should be purged with an inert gas. PF-06869206 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PF-06869206 in ethanol is approximately 10 mg/ml and approximately 1 mg/ml in DMSO and DMF.

PF-06869206 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-06869206 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. PF-06869206 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

PF-06869206 is an orally bioavailable inhibitor of sodium-phosphate cotransporter 2a (NaPi2a; IC<sub>50</sub> = 380 nM).<sup>1</sup> It is selective for NaPi2a over NaPi2b, NaPi2c, PiT-1, and PiT-2 (IC<sub>50</sub>s = >25 μM).

## Reference

1. Filipski, K.J., Sammons, M.F., Bhattacharya, S.K., *et al.* Discovery of orally bioavailable selective inhibitors of the sodium-phosphate cotransporter NaPi2a (SLC34A1). *Med. Chem. Lett.* **9**(5), 440-445 (2018).

### 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, 10/14/2022

## 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](http://WWW.CAYMANCHEM.COM)