

# PRODUCT INFORMATION

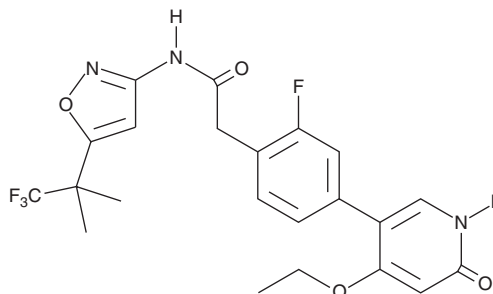


**GSK3179106**

Item No. 24023

**CAS Registry No.:** 1627856-64-7  
**Formal Name:** 4-(4-ethoxy-1,6-dihydro-6-oxo-3-pyridinyl)-2-fluoro-N-[5-(2,2,2-trifluoro-1,1-dimethylethyl)-3-isoxazolyl]-benzeneacetamide  
**Synonym:** RET Kinase Inhibitor 1

**MF:** C<sub>22</sub>H<sub>21</sub>F<sub>4</sub>N<sub>3</sub>O<sub>4</sub>  
**FW:** 467.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 233 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

GSK3179106 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK3179106 in the solvent of choice, which should be purged with an inert gas. GSK3179106 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK3179106 in ethanol is approximately 25 mg/ml and approximately 50 mg/ml in DMSO and DMF.

GSK3179106 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK3179106 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK3179106 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

GSK3179106 is an inhibitor of rearranged during transfection (RET) kinase (IC<sub>50</sub> = 0.4 nM).<sup>1</sup> It is selective for RET kinase over VEGF receptor 2 (VEGFR2/KDR; IC<sub>50</sub> = 109 nM). GSK3179106 reduces the number of abdominal contractions induced by colorectal distension in acetic acid-sensitized rats when administered at a dose of 10 mg/kg twice per day.

## References

1. Schenck Eidam, H., Russell, J., Raha, K., *et al.* Discovery of a first-in-class gut-restricted RET kinase inhibitor as a clinical candidate for the treatment of IBS. *ACS Med. Chem. Lett.* **9**(7), 623-628 (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, 12/22/2022

## 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)