

# PRODUCT INFORMATION



**BI-1935**

Item No. 41327

**CAS Registry No.:** 940954-41-6  
**Formal Name:** 1-(2-ethoxyethyl)-1,6-dihydro-6-oxo-N-[5-[3-(3-pyridinyl)-5-(trifluoromethyl)-1H-pyrazol-1-yl]-2-pyridinyl]-3-pyridinecarboxamide

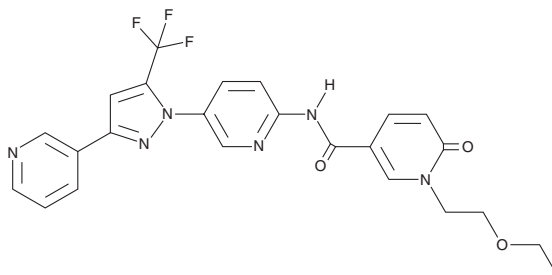
**MF:** C<sub>24</sub>H<sub>21</sub>F<sub>3</sub>N<sub>6</sub>O<sub>3</sub>  
**FW:** 498.5

**Purity:** ≥90%

**Supplied as:** A 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

BI-1935 is supplied as a solid. A stock solution may be made by dissolving the BI-1935 in the solvent of choice, which should be purged with an inert gas. BI-1935 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

BI-1935 is an inhibitor of soluble epoxide hydrolase (sEH; IC<sub>50</sub> = 7 nM). It is selective for sEH over the cytochrome P450 (CYP) isoforms CYP2J2, CYP2C9, and CYP2C19 (IC<sub>50</sub>s = >1,000 nM for all). See the Structural Genomics Consortium (SGC) website for more information.

### 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, 01/30/2026

## 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