

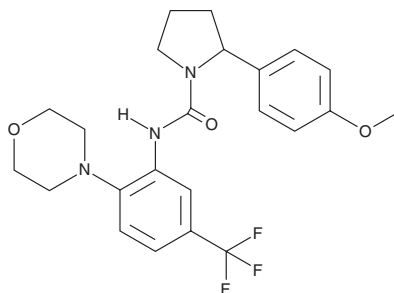
# PRODUCT INFORMATION



**ZINC40099027**

Item No. 42956

**CAS Registry No.:** 1211825-25-0  
**Formal Name:** 2-(4-methoxyphenyl)-N-[2-(4-morpholinyl)-5-(trifluoromethyl)phenyl]-1-pyrrolidinecarboxamide  
**Synonym:** ZN27  
**MF:** C<sub>23</sub>H<sub>26</sub>F<sub>3</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 449.5  
**Purity:** ≥98%  
**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

ZINC40099027 is supplied as a solid. A stock solution may be made by dissolving the ZINC40099027 in the solvent of choice, which should be purged with an inert gas. ZINC40099027 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

ZINC40099027 is an activator of focal adhesion kinase (FAK).<sup>1</sup> It induces FAK autophosphorylation in Caco-2 colon adenocarcinoma cells when used at a concentration of 10 nM. ZINC40099027 (10 nM) increases migration in a wound healing assay using Caco-2 cells. It increases wound closure and decreases ulcer area in a mouse model of acid-induced intestinal ulceration.

## Reference

1. Wang, Q., More, S.K., Vomhof-DeKrey, E.E., *et al.* Small molecule FAK activator promotes human intestinal epithelial monolayer wound closure and mouse ulcer healing. *Sci. Rep.* **9**(1), 14669 (2019).

### 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, 03/24/2025

## 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)