

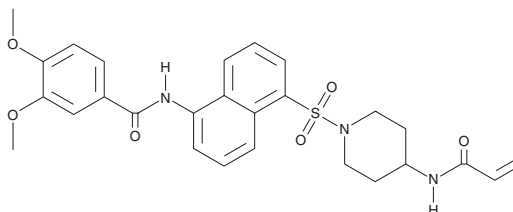
PRODUCT INFORMATION



BJJF078

Item No. 42397

CAS Registry No.: 2531244-56-9
Formal Name: 3,4-dimethoxy-N-[5-[[4-[(1-oxo-2-propen-1-yl)amino]-1-piperidinyl]sulfonyl]-1-naphthalenyl]-benzamide
MF: C₂₇H₂₉N₃O₆S
FW: 523.6
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

BJJF078 is supplied as a solid. A stock solution may be made by dissolving the BJJF078 in the solvent of choice, which should be purged with an inert gas. BJJF078 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

BJJF078 is an inhibitor of transglutaminase 2 (TG2; IC₅₀s = 41 and 54 nM for the human and mouse enzymes, respectively).¹ It is selective for TG2 over Factor XII (IC₅₀ = 22 μM for the human enzyme) but does not inhibit TG1 (IC₅₀ = 0.16 μM for the human enzyme).

Reference

1. Chrobok, N.L., Bol, J.G.J.M., Jongenelen, C.A., *et al.* Characterization of transglutaminase 2 activity inhibitors in monocytes *in vitro* and their effect in a mouse model for multiple sclerosis. *PLoS One* **13**(4), e0196433 (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, 12/16/2024

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