

PRODUCT INFORMATION



JB1-1044

Item No. 45845

CAS Registry No.: 2304852-21-7

Formal Name: [(3R)-3-amino-1-piperidinyl]
[2-[7-chloro-1-(cyclopropylmethyl)-1H-
indol-2-yl]-4,5-dihydroimidazo[1,5,4-
de][1,4]benzoxazin-8-yl]-methanone

MF: C₂₇H₂₈ClN₅O₂

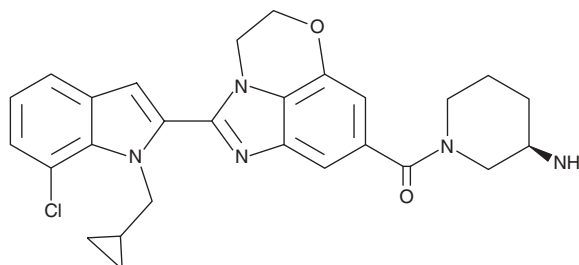
FW: 490.0

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

JB1-1044 is supplied as a solid. A stock solution may be made by dissolving the JB1-1044 in the solvent of choice, which should be purged with an inert gas. JB1-1044 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

JB1-1044 is an inhibitor of peptidyl arginine deiminase 4 (PAD4; IC₅₀ = 102 nM).¹

Reference

1. Luo, H., Zuo, H., Zhu, J., *et al.* Computer-aided discovery of peptidyl arginine deiminase 4 inhibitors to suppress NETosis and accelerate diabetic wound healing. *Chinese Chem. Lett.* **37**(6), 111510 (2026).

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, 07/13/2026

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