

PRODUCT INFORMATION

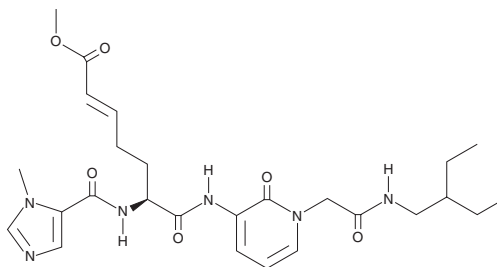


ZED-1227

Item No. 42333

CAS Registry No.: 1542132-88-6
Formal Name: (2E,6S)-7-[[1-[2-[(2-ethylbutyl)amino]-2-oxoethyl]-1,2-dihydro-2-oxo-3-pyridinyl]amino]-6-[[[(1-methyl-1H-imidazol-5-yl)carbonyl]amino]-7-oxo-2-heptenoic acid, methyl ester

Synonym: TAK-227
MF: $C_{26}H_{36}N_6O_6$
FW: 528.6
Purity: $\geq 95\%$
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

ZED-1227 is supplied as a solid. A stock solution may be made by dissolving the ZED-1227 in the solvent of choice, which should be purged with an inert gas. ZED-1227 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

ZED-1227 is an inhibitor of transglutaminase 2 (TG2; $IC_{50} = 0.053 \mu\text{M}$).¹ It is selective for TG2 over TG1, TG3, TG6, and Factor XIIIa₂ (IC_{50} s = 24.86, >50, 6.44, and >50 μM , respectively).

Reference

1. Büchold, C., Hils, M., Gerlach, U., *et al.* Features of ZED1227: The first-in-class tissue transglutaminase inhibitor undergoing clinical evaluation for the treatment of celiac disease. *Cells* **11**(10), 1667 (2022).

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, 12/04/2024

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