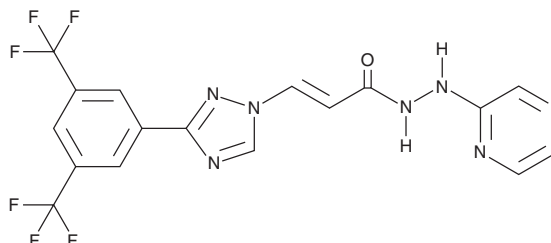


PRODUCT INFORMATION



(E)-KPT-330
Item No. 39562

CAS Registry No.: 1421923-86-5
Formal Name: 3-[3-[3,5-bis(trifluoromethyl)phenyl]-1H-1,2,4-triazol-1-yl]-2-(2-pyrazinyl)hydrazide, (2E)-2-propenoic acid
Synonyms: KTP-330, *trans*-Selinexor
MF: C₁₇H₁₁F₆N₇O
FW: 443.3
Purity: ≥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

(E)-KPT-330 is supplied as a solid. A stock solution may be made by dissolving the (E)-KPT-330 in the solvent of choice, which should be purged with an inert gas. (E)-KPT-330 is soluble in DMSO.

Description

(E)-KPT-330 is a metabolite of the exportin 1 (XPO1/CRM1) inhibitor selinexor.¹ It is formed from selinexor by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Mancuso, J. Selinexor (Xpovio), an XPO1 inhibitor and a new class of therapeutics for treating multiple myeloma. *Current Drug Synthesis*. Li, J.J., editor, 1st edition, John Wiley & Sons (2023).

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, 10/31/2023

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