

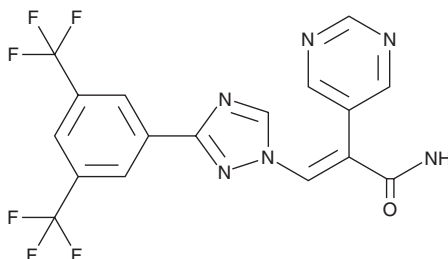
PRODUCT INFORMATION



Eltanexor

Item No. 43585

CAS Registry No.: 1642300-52-4
Formal Name: α E-[[3-[3,5-bis(trifluoromethyl)phenyl]-1H-1,2,4-triazol-1-yl]methylene]-5-pyrimidineacetamide
Synonym: KPT-8602
MF: C₁₇H₁₀F₆N₆O
FW: 428.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Description

Eltanexor is an inhibitor of exportin 1 (XPO1/CRM1).¹ It inhibits the binding of XPO1/CRM1 to various nuclear export signals (NESs) in cell-free pull-down assays when used at a concentration of 10 μ M. Eltanexor inhibits Cas9-encoding mRNA nuclear export in HEK293T cells in a concentration-dependent manner and CRISPR-Cas9 gene editing in a reporter assay using HEK293 cells (IC₅₀ = 94 nM).² It inhibits human cytomegalovirus (CMV) replication in primary human foreskin fibroblasts (IC₅₀ = 37.62 nM).³ Eltanexor (100 nM) induces I κ B α and NF- κ B p65 nuclear accumulation and inhibits RANKL-induced osteoclast differentiation in isolated mouse bone marrow-derived macrophages (BMDMs).⁴ It induces apoptosis in primary human chronic lymphocytic leukemia (CLL) cells and increases survival and decreases spleen leukemic burden in an MV4-11 acute myeloid leukemia (AML) mouse xenograft model.¹ Eltanexor (5 mg/kg five times per week) also increases locomotor activity and tibialis anterior myofiber size and decreases serum osteopontin (OPN) levels in the D2-*mdx* mouse model of Duchenne muscular dystrophy (DMD).⁵

References

1. Hing, Z.A., Fung, H.Y., Ranganathan, P., *et al.* Next-generation XPO1 inhibitor shows improved efficacy and *in vivo* tolerability in hematological malignancies. *Leukemia* **30**(12), 2364-2372 (2016).
2. Cui, Y.R., Wang, S.J., Ma, T., *et al.* KPT330 improves Cas9 precision genome- and base-editing by selectively regulating mRNA nuclear export. *Commun. Biol.* **5**(1), 237 (2022).
3. Liao, Y., Ke, X., Deng, T., *et al.* The second-generation XPO1 inhibitor eltanexor inhibits human cytomegalovirus (HCMV) replication and promotes type I interferon response. *Front. Microbiol.* **12**:675112, (2021).
4. Chen, J., Song, D., Xu, Y., *et al.* Anti-osteoclast effect of exportin-1 inhibitor eltanexor on osteoporosis depends on nuclear accumulation of I κ B α -NF- κ B p65 complex. *Front. Pharmacol.* **13**:896108, (2022).
5. English, K.G., Reid, A.L., Samani, A., *et al.* Next-generation SINE compound KPT-8602 ameliorates dystrophic pathology in zebrafish and mouse models of DMD. *Biomedicines* **10**(10), 2400 (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, 08/15/2025

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