

PRODUCT INFORMATION



GMX1777

Item No. 38798

CAS Registry No.: 432037-57-5
Formal Name: 4-[[[6-(4-chlorophenoxy)hexyl]imino](cyanoamino)methyl]amino]-1-(3-oxo-2,4,7,10,13,16-hexaoxaheptadec-1-yl)-pyridinium, monochloride

Synonyms: EB1627, Teglalinad

MF: C₃₀H₄₃ClN₅O₈ • Cl

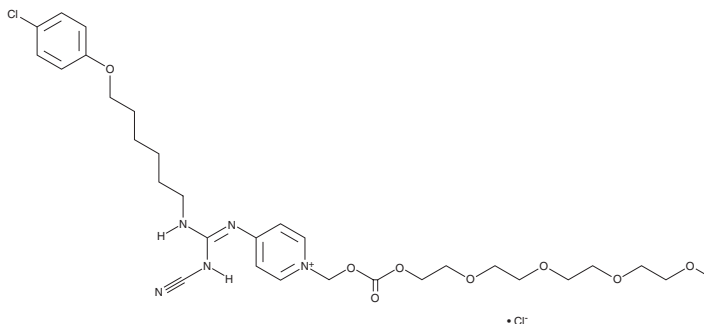
FW: 672.6

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

GMX1777 is supplied as a solid. A stock solution may be made by dissolving the GMX1777 in the solvent of choice, which should be purged with an inert gas. GMX1777 is soluble in DMSO.

Description

GMX1777 is a prodrug form of the nicotinamide phosphoribosyltransferase (Nampt) inhibitor CHS-828 (Item No. 11021).¹ Intravenous administration of GMX1777 (75 and 150 mg/kg) reduces NAD⁺ levels in tumors, but not muscle tissue, in a DMS 273 lung cancer mouse xenograft model. GMX1777 (50 mg/kg per day), in combination with radiation, decreases tumor diameter and microvessel density in C666-1 nasopharyngeal and FaDu head and neck squamous cell carcinoma (HNSCC) mouse xenograft models.²

References

1. Beuparlant, P., Bédard, D., Bernier, C., *et al.* Preclinical development of the nicotinamide phosphoribosyl transferase inhibitor prodrug GMX1777. *Anticancer Drugs* **20**(5), 346-354 (2009).
2. Kato, H., Ito, E., Alajez, N.M., *et al.* Efficacy of combining GMX1777 with radiation therapy for human head and neck carcinoma. *Clin. Cancer Res.* **16**(3), 898-911 (2010).

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, 05/26/2023

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