

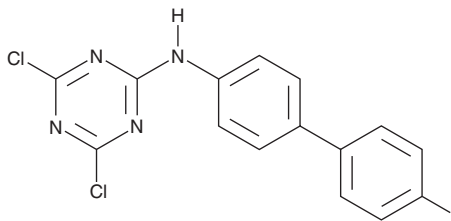
PRODUCT INFORMATION



KEA1-97

Item No. 42049

CAS Registry No.: 2138882-71-8
Formal Name: 4,6-dichloro-N-(4'-fluoro[1,1'-biphenyl]-4-yl)-1,3,5-triazin-2-amine
MF: C₁₅H₉Cl₂FN₄
FW: 335.2
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

KEA1-97 is supplied as a solid. A stock solution may be made by dissolving the KEA1-97 in the solvent of choice, which should be purged with an inert gas. KEA1-97 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

KEA1-97 is a dichlorotriazine (DCT) inhibitor of the protein-protein interaction between thioredoxin (Trx) and caspase-3.¹ It binds to lysine 72 (K72) of Trx and inhibits DCT-alkyne labeling of recombinant Trx at K72 (IC₅₀ = 10 μM). KEA1-97 (100 μM) inhibits the interaction between Trx and caspase-3 in a cell-free assay. It reduces survival and proliferation of and induces apoptosis in Trx-overexpressing 231MFP breast cancer cells. *In vivo*, KEA1-97 (5 mg/kg) reduces tumor volume in a 231MFP mouse xenograft model.

Reference

1. Anderson, K.E., To, M., Olzmann, J.A., *et al.* Chemoproteomics-enabled covalent ligand screening reveals a thioredoxin-caspase 3 interaction disruptor that impairs breast cancer pathogenicity. *ACS Chem. Biol.* **12**(10), 2522-2528 (2017).

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, 11/08/2024

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