PRODUCT INFORMATION



Lys05 (hydrochloride)

Item No. 27334

CAS Registry No.:	1391426-24-6	
Formal Name:	N ² -(7-chloro-4-quinolinyl)-N ¹ -[2-[(7-chloro-	
	4-quinolinyl)amino]ethyl]-N ¹ -methyl-1,2- ethanediamine, trihydrochloride	H_N_N_H
Synonym:	Lys01	Ĩ
MF:	$C_{23}H_{23}CI_2N_5 \bullet 3HCI$	
FW:	549.8	
Purity:	≥98%	
UV/Vis.:	λ _{max} : 221, 257, 292, 331, 344 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	• 3001
Stability:	≥4 years	

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

Laboratory Procedures

Lys05 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the lys05 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Lys05 (hydrochloride) has an approximate solubility of 2 mg/ml in DMSO. Lys05 (hydrochloride) is slightly soluble in ethanol.

Lys05 (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Lys05 is an inhibitor of autophagy.¹ It induces p62 accumulation and increases the ratio of LC3-II to LC3-I in C8161 melanoma cells in a concentration-dependent manner. Lys05 (76 mg/kg) increases the number of autophagic vesicles per cell in tumors in a C8161 mouse xenograft model and reduces tumor volume and growth in a 1205Lu melanoma mouse xenograft model. It also reduces tumor growth rate, volume, and weight in an HT-29 colon cancer mouse xenograft model when administered at doses greater than or equal to 10 mg/kg but increases bowel thickness and obstruction when administered at 80 mg/kg.

Reference

1. McAfee, Q., Zhang, Q., Samanta, A., et al. Autophagy inhibitor Lys05 has single-agent antitumor activity and reproduces the phenotype of a genetic autophagy deficiency. Proc. Natl. Acad. Sci. U.S.A. 109(21), 8253-8258 (2012).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 12/15/2022

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